

Discovery Life Sciences
6 Henshaw Street
Woburn, MA 01801
Tel: (866) 838-2798
info@dls.com
<https://www.dls.com/>

UltraPool™ HLM 150

Catalog No.	452117
Lot No.	38298
Qty./Package	0.5 mL
Protein Content	20 mg/mL in 250 mM sucrose
Storage Conditions	Store at -80°C
Release Date	2022 October
Expiration Date	2032 September

Assay Results

Enzyme Measured	Assay	Enzyme Activity [in pmol/(mg x min)]
Total P450	Omura and Sato	350 pmol/mg
OR	Cytochrome c Reductase	300 nmol/(mg x min)
Cytochrome b ₅	Spectrophotometric	400 pmol/mg
CYP1A2	Phenacetin O-deethylase	460
CYP2A6	Coumarin 7-hydroxylase	700
CYP2B6	Bupropion hydroxylase	300
CYP2C8	Paclitaxel 6 α -hydroxylase	150
CYP2C9	Diclofenac 4'-hydroxylase	2700
CYP2C19	(S)-Mephenytoin 4'-hydroxylase	45
CYP2D6	Bufuralol 1'-hydroxylase (The amount of activity inhibited by 1 μ M quinidine.)	110
CYP2E1	Chlorzoxazone 6-hydroxylase	2100
CYP3A4	Testosterone 6 β -hydroxylase	4100
CYP4A11	Lauric acid 12-hydroxylase	1700
FMO	Methyl p-Tolyl Sulfide Oxidase	700
UGT1A1	Estradiol 3-Glucuronidation	1100
UGT1A4	Trifluoperazine N- Glucuronidation	620
UGT1A6	Serotonin Glucuronidation	21000
UGT1A9	Propofol Glucuronidation	1800
UGT2B7	AZT Glucuronidation	1600
Western Blot Results		
CYP3A4	Western Blot	91 pmol/mg
CYP3A5	Western Blot	40 pmol/mg

- All cytochrome P450 assays were conducted at 0.8 mg/mL protein (except CYP2B6 which was at 0.1 mg/mL, CYP2C19 which was at 0.3 mg/mL, and CYP3A4 which was at 0.5 mg/mL) with a NADPH generating system (1.3 mM NADP, 3.3 mM glucose 6-phosphate, 3.3 mM MgCl₂, and 0.4 U/mL glucose 6-phosphate dehydrogenase) and incubated for 20 minutes (except CYP2B6 which was incubated for 5 minutes and CYP2C8, CYP2C9, CYP2C19, CYP3A4 and CYP4A11 which were incubated for 10 minutes). 0.1 M Potassium phosphate buffer, pH 7.4 was used for all P450 enzymes (except 0.5 M Potassium phosphate buffer, pH 7.4 was used for CYP2B6 and CYP2C19 and 0.1 M Tris buffer, pH 7.5 was used for CYP2A6, CYP2C9 and CYP4A11). The FMO assay was conducted at 0.8 mg/mL protein with the same NADPH generating system, 1.2 mM diethylenetriaminepentacetic acid, and 0.5 mg/mL Triton X-100 in 0.05 M glycine buffer, pH 9.5, and was incubated for 10 minutes. UGT Glucuronidation assays were conducted at 0.5 mg/mL protein for UGT1A1 and UGT1A4, 0.1 mg/mL for UGT1A6, 0.15 mg/mL for UGT1A9, and 0.8 mg/mL for UGT2B7 with 2 mM UDPGA, 10 mM MgCl₂, and 25 μ g/mL Alamethicin in 0.1 M Tris buffer, pH 7.5. UGT1A1 and were incubated for 30 minutes, UGT1A4 for 20 minutes, UGT1A6 for 15 minutes, UGT1A9 for 10 minutes, and UGT2B7 for 25 minutes. Activities expressed as pmol product per mg protein x minute (except cytochrome c reductase which is expressed as nmol product per mg protein x minute).
- The Western Blot assay was carried out using standard protocols. SDS-gel electrophoresis was by the method of Laemmli (Laemmli, U.K, 1970, *Nature*, 227: 680-685.). CYP protein abundance in HLM was quantitated using authentic standards derived from P450 isoforms.
- The pool is comprised of equal milligrams of microsome per donor.
- The protein concentration reported has been demonstrated to be within an acceptable range of 18-24 mg/mL. Range reported is due to intrinsic variability in the assay.

For Research Use Only. Not for use in diagnostic or therapeutic procedures.

© Discovery Life Sciences

Discovery Life Sciences
6 Henshaw Street
Woburn, MA 01801
Tel: (866) 838-2798
info@dls.com
<https://www.dls.com/>

UltraPool™ HLM 150

Donor Information

Specimen	Gender	Age (years)	Race	Cause of Death	Social History	Medical History	Medication given during Hospitalization
*HMC1003	Male	32	C	Other 2 nd Cardiovascular	Cigarettes: more than 1 PPD x 15 yrs; Quit for 1 yr 5-6 yrs ago. ETOH: (Beer) 1-2 beers a wk since 21 yrs old. Marijuana, Sleep pills: Smoked unknown amounts for an unknown duration and frequency.	Heart Disease; Unnamed diagnosis on admission.	Ativan, Heparin, and Ancef. Flushed and stored in SPS-1.
HMC1004	Male	39	C	Anoxia 2 nd Drug Intoxication	Cigarettes: 1.5 PPD since 18 yrs old. ETOH: (Distilled Spirits) Drank unknown amount for unknown duration. Heroin: Drug was taken by any method possible; Unknown amount and frequency for 3 months.	HTN: Diagnosed ~ 2 yrs ago; Compliant w/ treatment. Depression: Diagnosed with clinical depression and anxiety; unknown when diagnosed or if compliant with treatment. Prilosec and other medication were taken at home, but they were not named. (as perNOK).	Heparin, Regitine, Versed, Fentanyl, Solumedrol, Furosemide and Insulin. Flushed and stored in UW solution.
HFC1005	Female	49	C	Anoxia 2 nd Natural Causes	Cigarettes- 5 cigarettes per day X 25 yrs; Quit in 2009. ETOH (Beer, Wine) 1-2 drinks per month for an unknown duration.	Pulmonary fibrosis; Diagnosed during hospital stay; Unknown if compliant w/ treatment. No meds taken at home (as per NOK). Pneumonia; diagnosed a yr prior and never really recovered.	Insulin, Decadron, and Heparin. Flushed and stored in SPS-1.
*HMC1007	Male	61	C	HT 2 nd Natural Causes	ETOH (Beer): 3-4 drinks per wk x 40 yrs.	Asthma: Childhood asthma; No symptoms or treatment needed in yrs. No meds take at home (as per NOK).	Heparin, KCL, and MgSO4. Flushed and stored in SPS-1.
HFC1009	Female	24	C	HT 2 nd Blunt Injury	ETOH (Wine) 3 wine coolers at a time, 3 X during the summer since age 21.	No meds taken at home (as per NOK).	Levophed, Albuterol, Lasix, Insulin, Antihypertensives, Heparin, CaGluconate, Lopressor, Cefazolin, Flagyl, Vancomycin, Cefepime, Duoneb, Keppra, Mannitol, KPhos, NaPhos, Heparin. Flushed and stored in HTK.
*HFC1013	Female	40	C	CVA 2 nd ICH	ETOH; Social drinker; Unknown type, amount, frequency, or duration.	MVI (as per NOK). Pt had no previous medical history aside from child birth.	Levophed, Neosynephrine, Hydrocortisone, NaBicarb, Insulin, NaCl, NaHCO3, Solumedrol, DDAVP, Zosyn. Flushed and stored in UW Solution.
*HFC1017	Female	63	C	Other 2 nd ICH	Non-smoker; Non-drinker	The donor did not take any OTC or prescribed medication on a regular basis (as per NOK).	Levophed, Solumedrol, Ancef, Hespan, PRBCs. Flushed and stored in SPS-1.

For Research Use Only. Not for use in diagnostic or therapeutic procedures.

© Discovery Life Sciences

Discovery Life Sciences
 6 Henshaw Street
 Woburn, MA 01801
 Tel: (866) 838-2798
info@dls.com
<https://www.dls.com/>

UltraPool™ HLM 150

Donor Information

Specimen	Gender	Age (years)	Race	Cause of Death	Social History	Medical History	Medication given during Hospitalization
HMC1020	Male	50	C	Anoxia 2 nd Cardiovascular	Cigarettes: 6 cigarettes PPD x 30 yrs on and off; Chewing Tobacco 30 yrs ago; Used for 2 yrs and the quit. ETOH (Beer) During his 20's, 12 beers / weekend; During the last few yrs, 6 beers / month since age 10. Marijuana; (Inhaled) Unknown amount used weekly as a teenager; Crystal meth (Inhaled) 3-4 x / week; Unknown amount; Used from age 25-50; Was currently using; Cocaine (Intranasal) 1-2 x / month; Unknown amount; Used from age 20-50; Last ime used was 25.	Asthma; Diagnosed as a child: Resolved as an adult. DM x 15 yrs; Unknown type; Non-compliant with meds except while incarcerated. No meds taken at home (as per NOK)	Vancomycin, Vasopressin, Zosyn, NaHCO ₃ , Propofol, Levophed, Insulin, Albuterol, NS, Solumedrol, Kayexalate, Hespan, CaCl, Heparin, Mannitol, Lasix. Flushed and stored in UW solution.
*HFC1021	Female	46	C	Pneumonia 2 nd Natural Causes	ETOH (Wine, Wine Coolers) 1-2 drinks during special occasions; Unknown duration.	IBS; It was cleared up once Gall Bladder was removed; Bloody Stools occasionally-last episode was a couple of months ago. Frequent Bladder Infections over 20 yrs. Ago. T4, Premarin, Glucosamine (as per NOK).	Kayexalate, Acetaminophen, KCl, Amiodarone, Acetazolamide, Lasix, MgSulfate, Lovenox, Insulin, TPN, T4, Rocephin, Levaquin, NaBicarb, Hydrocortisone, Vancomycin, Levophed, Nimbex, Cisatracurium, Propofol, Merrem, Solumedrol. Flushed and stored in SPS-1.
*HMC1023	Male	48	C	Anoxia	Cigarettes: 1/4 PPD x 30 yrs. ETOH: 2 cases of Beer per week, unknown duration. Cocaine (Inhaled, Intranasal) 20 yrs ago, Unknown amount, duration, or frequency; Marijuana (Smoked) currently, Unknown amount, duration, or frequency.	HTN x 2 yrs; Compliant w/ meds. Liprosil, Men's daily vitamin (as per NOK).	Levophed, Solumedrol, Epinephrine, Vasopressin, Insulin, Dilantin, Magnesium Sulfate, Potassium Chloride, Dextrose, Ciprofloxacin, Vancomycin, Zosyn, Methylprednisone, Levothyroxine. Flushed and stored in SPS-1.

For Research Use Only. Not for use in diagnostic or therapeutic procedures.

© Discovery Life Sciences

Discovery Life Sciences
6 Henshaw Street
Woburn, MA 01801
Tel: (866) 838-2798
info@dls.com
<https://www.dls.com/>

UltraPool™ HLM 150

Donor Information

Specimen	Gender	Age (years)	Race	Cause of Death	Social History	Medical History	Medication given during Hospitalization
HMC1027	Male	66	C	Head Trauma 2 nd ICH	Cigarettes: 2.5 PPD X 35 yrs; quit in 2003. ETOH: occasional beer, but has not done in yrs.	Coronary Artery Disease; stents x 4; believed to be due to exposure to Agent Orange in Vietnam. Bipolar disorder, Post Traumatic Stress Disorder, Dementia x 2-3 yrs and Confusion/memory loss due to Dementia. Yellow Jaundice due to Lithium toxicity 6 yrs ago that caused acute liver failure. Failure was resolved; unsure of treatment. Acute Renal Failure 6 yrs ago due to Lithium Toxicity. Acetaminophen, Aspirin, Atorvastin, Benzotropine Mesylate, Donepezil, HCL, Olanzapine, Omeprazole, Caroxerine, Tamsoulin, and Vitamin D3 on a regular basis (as per NOK).	Levophed, Neosynephrine, Levothyroxine, T4, DDAVP, Hypertensives, Magnesium, Insulin, Ancef, Solumedrol, Lopressor, Cardizem and Platelets. Flushed and stored in HTK.
*HFC1028	Female	46	C	CVA 2 nd ICH	Cigarettes- 1 PPD X 28 yrs; Quit 3 months ago. ETOH- 1 drink of mixed liquor per week X 10 yrs; quit 7 yrs ago.	The donor suffered with spontaneous pneumothoraces X 2 in 2007; required surgical repair. Seizures and depression. Severe malnutrition due to Anorexia; recent PEG tube placement. Donor took Dilantin, Keppra, Levothyroxine, Folic Acid, Culturelle, Augmentin, Seroquel, and Oxycontin on a regular basis (as per NOK). Hypothyroidism and chronic migraines.	Levophed, Vasopressin, Dobutamine, Insulin, Vasopressin, Zosyn, Vancomycin, Albumin, Heparin and Phentolamine. Flushed and stored in HTK.
HMC1030	Male	32	C	Anoxia			Dopamine, T4, Insulin, Flushed and stored in HTK.
*HFC1035	Female	59	C	CVA 2 nd ICH	cigarettes- 0.5-1 PPD x 43 yrs. ETOH (Beer) 1 drink/ year X 30 yrs. Marijuana (Inhaled) 1 joint/ year X 30 yrs.	HTN; Diagnosed about 10 yrs ago, Compliant with treatment and medication. Intestinal issues (Diarrhea and Bloody Stools); Diagnosed within the last couple of yrs. Son reported it was getting better. Prednisone, Aspirin, Lopressor, Lisonpril, Pregabalin, Bupropin (meds taken at home as per NOK)	Esmolol, Solumedrol, Insulin, Antihypertensives, DDAVP, KCl, Labetalol, Nipride. Flushed and stored in SPS-1.
HMC1036	Male	56	C	HT 2 nd Blunt Injury	Cigarettes; 1 ppd X 40+ yrs; Chewing Tobacco, unknown amount X 4 yrs. ETOH (Beer) 3-4 daily X 40+ yrs. Crack (Inhaled) Unknown amount or frequency X 9 yrs; Marijuana (Inhaled) Unknown amount, frequency or duration.	As per donor's spouse donor did not have HTN; but as per donor's daughter he have HTN; Unknown when diagnosed or if compliant with meds. Unknown names of medications taken at home as reported by NOK.	Heparin, Fentanyl, Ativan, Levophed, Dextrose, DDAVP. Flushed and stored in SPS-1.

For Research Use Only. Not for use in diagnostic or therapeutic procedures.

© Discovery Life Sciences

Discovery Life Sciences
6 Henshaw Street
Woburn, MA 01801
Tel: (866) 838-2798
info@dls.com
<https://www.dls.com/>

UltraPool™ HLM 150

Donor Information

Specimen	Gender	Age (years)	Race	Cause of Death	Social History	Medical History	Medication given during Hospitalization
*HMC1040	Male	61	C	CVA 2 nd ICH	Cigarettes: 2 PPD x 25+ yrs. Marijuana (Inhaled) Unknown amount x 25 yrs, unknown when last used	No meds taken at home (as per NOK).	Dopamine, Neosynephrine, Vasopressin, Solumedrol, T4, Insulin, Antihypertensives, Cefazolin, Vancomycin, Zosyn, Heparin, Mannitol, Lasix, Cardene, Pitressin, Albumin, Keppra, Sodium BiCarbonate, LR, Ancef, KCl, KPhos, Flushed and stored in SPS-1.
*HFC1043	Female	53	C	Anoxia 2 nd Cardiovascular	Cigarettes; 0.5 PPD x 20 yrs. ETOH (Beer, Wine) Unknown amount, socially, unknown duration. Marijuana (Inhaled) Unknown amount, frequency, or duration.	HTN; Diagnosed 15-20 yrs ago, compliant with meds. Skin Cancer- Treatment included surgical removal 3 weeks ago. MVI, HTN, Pain and Anxiety (Meds taken at home as per NOK).	Heparin, Levophed, Vasopressin, Cefepime, Zosyn, Amiodarone. Flushed and stored in UW solution
*HFC1046	Female	57	C	CVA 2 nd ICH	ETOH (Scotch) 2-3 drinks/ week socially x 40 yrs.	Mitral Valve Prolapse; Unknown when diagnosed. Atenolol, MVI (meds taken at home per NOK).	Neosynephrine, Levophed, Solumedrol, Heparin, Ancef. Flushed and stored in UW solution
HMC1048	Male	36	C	CVA 2 nd ICH	ETOH (Beer, Whiskey) Social drinking 2-3 times per week-amount would fluctuate between 2-10 drinks per outing, since age 18. Cocaine - used since early 20's, NOT IV, use slowind down in the last 6 months, about 2 xs per month.	Stomach Ulcers on and off for the last year causing diarrhea. Steroid for Ulcers for 10 days, Metamucil (as per NOK).	Heparin, Mannitol, Rocephin, Azithromycin, Norepinephrine, Sodium Chloride, Epinephrine, Vasopressin, Dextrose, Solumedrol, T4, Insulin. Flushed and stored in SPS-1.
HMC1052	Male	26	C	Anoxia 2 nd Cardiovascular	Non-smoker; Non-drinker	Mental retardation. MVI, Zonegran, Trazadoen, Cogenten (as per NOK).	Versed, Kphos, Keppra, Vancomycin, Zosyn, Propofol, Heparin, Phenytoin, Levophed, Neosynephrine. Flushed and stored in UW solution.
HMC1055	Male	64	C	HT 2 nd Blunt Injury	ETOH (Beer, Wine) 1 beer a day, 5-7 days a week x 5 years. Occasional glass of wine.	HTN; Unknown when diagnosed; compliant with treatment. Lisinopril (as per NOK).	Heparin, Levophed, Neosynephrine, Dextrose, Solumedrol, Lasix, T4, Insulin, DDAVP. Flushed and stored in SPS-1.
*HMC1056	Male	56	C	CVA 2 nd ICH	Cigarettes: 1.5 PPD x 40 yrs. ETOH (Beer) 10 beers, three times a week x 40 yrs. Cocaine & Marijuana once a year, 10+ yrs ago.	Blockage in heart - Pacemaker, stent placed. Heart cath performed yearly. HTN; Diagnosed 10 yrs ago. DM; Diagnosed 10 yrs ago, treated with Diet. HTN meds, Lasix, Zocor, Beta pace, Coreg, Lisinopril, Inspra, ASA, and Eliquis (as per NOK).	Levophed, DDAVP, Rosephin, KCL. Flushed in UW, stored in HTK.
*HFC1058	Female	47	C	Anoxia 2 nd Natural Causes	Champagne; 1 glass every 6 weeks X 20 yrs.	Asthma; Unknown when diagnosed. HTN; Diagnosed 20+ yrs ago. DM Diabetes; Diagnosed 22 yrs ago. Gestational Diabetes in 1990. Treated by medication, Insulin, Lantis, Januvia.	Levophed, Vasopressin, Insulin, Propofol, Zosyn, Decadron, Heparin. Flushed and stored in SPS-1.

For Research Use Only. Not for use in diagnostic or therapeutic procedures.

© Discovery Life Sciences

Discovery Life Sciences
6 Henshaw Street
Woburn, MA 01801
Tel: (866) 838-2798
info@dls.com
<https://www.dls.com/>

UltraPool™ HLM 150

Donor Information

Specimen	Gender	Age (years)	Race	Cause of Death	Social History	Medical History	Medication given during Hospitalization
HMC1060	Male	55	C	HT 2 nd Blunt Injury	Cigarettes: 0.5 PPD x 23 yrs. Quit for 7 yrs then resumed smoking to present day. Chronic alcoholism. Up to a liter a Vodka a day. Has drank mouth wash when out of money & able to shop lift. Pt was in detox 3 weeks ago. Marijuana (Inhaled) unknown amount, daily x 50 yrs. Cocaine and Amphetamines casually. Last use of drugs was 10 yrs ago.	Chronic Bronchitis. Bipolar Disorder - stopped taking medication 6 months ago. Depression, unknown duration. Chronic Pancreatitis. Flare up 2 months ago. No meds at home per NOK in the DRAI.	Artificial tears, DuoNeb, Cipro, Pepcid, Calcium Gluconate, Heparin, Keppra, Insulin, Kphos, Magnesium Sulfate, Labetalol, Ativan, KCl, Hydralazine, Tylenol, Lasix, D5, Nitroprusside, Dextrose, Lasix, Antihypertensives, Albuterol. Flushed and stored in SPS-1.
HMC1062	Male	55	C	CVA 2 nd ICH	Cigars; 2 cigars, once a day x 3 yrs. ETOH (Beer, Liquor) 2 drinks, daily since age 14. Cocaine, THC - unknown amount for a year, last used 25 yrs ago.	HTN; Diagnosed 4 mos ago. Lisinopril, MVI (as per NOK).	Versed, Cardene, Tylenol, Vasotec, Neosynephrine, NS, D5W, Heparin. Flushed in HTK and stored in Viaspan.
HMC1066	Male	46	C	Anoxia 2 nd Drug Intoxication	Cocaine, Marijuana- unknown amount, regularly since age 13.	Fibromyalgia meds, MSRA+ meds, Dietary supplements (as per NOK).	Vasopressin, Lasix, NS, Zosyn, Protonix, Keppra, Potassium Chloride, DDAVP. Flushed and stored in SPS-1.
*HFC1067	Female	56	C	Anoxia 2 nd Drug Intoxication	Cigarettes- 1 PPD x 40 yr. Vodka; a pint twice a week since the age of 18.	For the past 2 yrs. Pt has been very forgetful. Albuterol, Xanax, Azithromycin, Benzonatate, Citalopram, Lamictal, Tramadol, Trazadone, Zyrtec (as per NOK).	Versed, Morphine, D5W, Epinephrine, BiCarb, Levophed, Dopamine, Insulin, Zosyn, Solumedrol, Vancomycin, T4, D50, Vitamin K, Magnesium Sulfate, Calcium Gluconate, Albumin. Flushed and stored in SPL-1.
HMC1070	Male	55	C	Anoxia 2 nd Cardiovascular	Beer: one can, once a month for the past 5 yrs.	Asthma as a child, but cleared as he got older. He now has pulmonary fibrosis since 2014. Esbriet, Naprosyn, Zolofit until June 20115, MVI, Glucosamine (as per NOK).	Levophed, Albuterol, DuoNeb, Precedex, Prednisone, Solumedrol. Lasix, Bumex. Flushed and stored in HTK.
*HMC1071	Male	50	C	Anoxia 2 nd Drowning	Cigars; 3 a yr x 20 yrs. (Beer) 6-12 beers a day x 30 yrs.	Normal healthy	Heparin, Morphine, Cardene, Labetalol, Hydralazine, Insulin, Rocephin, Metoprolol, Phenytoin, Albuterol. Flushed and stored in SPS-1.
*HMC1073	Male	72	C	Anoxia	Cigarettes: 1 PPD x 20 yrs. ETOH: couple cans of beer on the weekends since 20 yrs old. Quit drinking 15 yrs ago.	DRAI, HTN; unknown duration, Asthma; unknown onset or duration, Kidney Disease; unknown duration, details or treatment, IDDM > 10 yrs., carpal tunnel, sleep apnea. HTN meds taken at home per NOK. Compliant w/ tx.	Levophed. Flushed and stored in SPS-1.

For Research Use Only. Not for use in diagnostic or therapeutic procedures.

© Discovery Life Sciences

Discovery Life Sciences
 6 Henshaw Street
 Woburn, MA 01801
 Tel: (866) 838-2798
info@dls.com
<https://www.dls.com/>

UltraPool™ HLM 150

Donor Information

Specimen	Gender	Age (years)	Race	Cause of Death	Social History	Medical History	Medication given during Hospitalization
*HMC1076	Male	61	C	ICH-Stroke		HTN > 10 yrs. PMHX: CVA- 8 months ago-untreated, high blood pressure, Type II DM, Cholecystectomy, CAD, high cholesterol. Blood pressure meds; had not taken in past 2 yrs due to no insurance. Half a regular adult aspirin once a day. Gall bladder surger 25 yrs ago. Previous stroke. On HCTZ for HTN; non-compliant with meds for past 3 yrs due to no insurance.	Heparin. Flushed and stored in HTK.
HMC1078	Male	65	C	ICH-Stroke	Cigarettes: 1 PPD x 35 yrs. ETOH: alcohol on weekends; 4 beers x 44 yrs.	Normal healthy	Levophed, Neosynephrine, Solumedrol, Lasix, Duonebs, Rocephin, Protonix, 0.45NS, CaCl, NaHCO3,
*HMC1080	Male	56	C	ICH-Stroke	Cigarettes: 2 PPD x 15 yrs; Quit 15 yrs ago. Drank socially on Christmas.	High Blood pressure and high cholesterol; Diagnosed 10 yrs ago. HTN > 10 yrs.; treatment compliant, HDL. Vertigo x 1 wk with headache, hearing loss in both ears, and tinnitus. Kidney Stones 20 yrs ago; no problems since. Benzapryl, hydrochlorothorizide, norvasc, coreg, and lotensin.	Atropine, Adensoine, Heparin, Arganine vasopressin,
*HFC191	Female	65	C	Anoxia	Tobacco use: 1.5 PPD x 20 yrs, quit 15 yrs ago.	Asthma, Emphysema and COPD for the last 15 yrs. Cardiac disease; diagnosed 15 yrs prior. Diabetes; diagnosed 5 yrs prior; diet regulated for 3 yrs, on meds last 2.	Protonix, Levaquin, Rocephin, Regular Insulin, Albumin, Vasopressin, Ancef, Zosin, D50, Solumedrol, Levothyroxin, Levo drip and Narcan
*HMC199	Male	42	C	Head Trauma	Cigarettes; 1/2 PPD x 23 yrs. 1-2 drinks q wk since age 21. Marijuana; one joint approx 6 times q yr. Family denied Cocaine use even though pt tested positive on admission.	Colon Sphincter repair 5-6 yrs ago. Large liver with 5% - 10% fat content by visual inspection.	Dilantin, Insulin, Kcl, Mannitol, Vancomysin, Etomidate, Lidocaine, Tetanus, Vecuronium.
HMC254	Male	24	C	Anoxia	Cigarettes; 1 1/2 PPD since 16. Beer; 12 pck over the weekend a 2-3 wks. Meth; approx 6-7 yrs ago x 3-4 yrs; amount and frequency unknown. Marijuana; every 3-6 months. Painkillers; Lortab (when he could get it), donor did not do any IV drugs.	Cardiac Ablasion d/t rapid heart beat (similarly to panic attack) 3-4 yrs ago. Enlarged Spleen and over production of blood. Xanax, Prozac and Thioridazine (as per NOK).	Dopamine, Neosynephrinem and Solucortef. Narcan and Romazicon administered

For Research Use Only. Not for use in diagnostic or therapeutic procedures.

© Discovery Life Sciences

Discovery Life Sciences
6 Henshaw Street
Woburn, MA 01801
Tel: (866) 838-2798
info@dls.com
<https://www.dls.com/>

UltraPool™ HLM 150

Donor Information

Specimen	Gender	Age (years)	Race	Cause of Death	Social History	Medical History	Medication given during Hospitalization
HMC275	Male	54	C	CVA	Cigarettes- 2 PPD X 20 yrs. Rum and Coke- nightly and Wine coolers > 25 yrs ago; no regular drinking with in 15 yrs.	Asthma- treated with inhaler; diagnosed 5 yrs ago. Hypertension; Diagnosed 5 yrs ago; Compliant w/ meds. Diabetes; Diagnosed 5 yrs ago; non-insulin dependent. Pt was treated for Hepatitis in his late teens-type in unk. Blood was found in pt's urine in Feb. 2008. HTN, Cholesterol and Diabetes meds (non insulin), asthma inhaler and nasal spray (as per NOK).	Morphine, Levophed, Neosyneprine, Dopamine, Vasopressin, Kphos, Ancef, solumedrol, and Cleocin, Flushed with UW solution.
*HFC289	Female	65	C	CHI	Cigarettes- 1-2 q d; duration unknown but pt quit since 1989. Wine- 1 glass (since adulthood on a rare basis.)	PT had cervical CA 20 yrs ago. leffexor (as reported by NOK.)	Insulin, Phospherytoin, DDAVP, Hydrocortisone, MGSo4, Mannitol, Sodium bicarb, Epinephrine, Fentanyl, Phenylephrine, Levothyroxine, Atomadate, SCCS (DA medication), Versed, Vecuronium, T4, Pitressin, NS, Zosyn and Vasopressin. Insitu: LR/UW and stored in UW solution.
*HMC290	Male	31	C	Head Trauma- Blunt Injury	Cigarettes- 1 PPD X 10 yrs. Beer/Liquor 2-3/ day X 4 yrs.		Levophed, D5w, 0.95 NaCL, FFP Plasma, and Solumedrol. Flushed and stored in HTK. Acinobactor, Ancef, and Clindamycin. Sedated on Propofol PRN and Fentanyl.
*HFC291	Female	29	C	Anoxia-GSW	Mixed Drinks- Rarely X 12 yrs. No tobacco	Gallbladder removed 2 yrs prior. C-section one yr prior.	Crystalloids, Insulin, Propofol, KCL, Levaquin, Lovenox, Levophed, 1/2 NS, and Magnesium Sulfate. Flushed and stored in HTK
HMC293	Male	55	C	Anoxia	Beer- Rarely.	Coronary Artery Disease. Heart Meds and Cholesterol Meds (as per NOK). History of previous MI and stent placement.	Insulin, Versed, Moxifloxacin, Albuterol, Keppra, D5NS, Cyrstalloids, Levophed, and Heparin. Epinephrine and atropine were administered. Flushed and stored in UW solution.
HFC300	Female	63	C	Anoxia	Beer- 3/day x 30 yrs.	Diabetes; Diagnosed 20yrs ago. Tx w/ Insulin Injections x 15 yrs.	Levophed, NS, Dopamine, Neosyneprine, Heparin, Insulin, Aztreonam, Decadron, Ancef, Nystatin, Erythromycin, Neomycin, Clindamycin, and Magnesium.

For Research Use Only. Not for use in diagnostic or therapeutic procedures.

© Discovery Life Sciences

Discovery Life Sciences
 6 Henshaw Street
 Woburn, MA 01801
 Tel: (866) 838-2798
info@dls.com
<https://www.dls.com/>

UltraPool™ HLM 150

Donor Information

Specimen	Gender	Age (years)	Race	Cause of Death	Social History	Medical History	Medication given during Hospitalization
HFC303	Female	59	C	ICH-Stroke	Non-smoker; Non-drinker	Cardizem (as per NOK)	Crystalloids, NS, Levophed, Albuterol, Furosemide, Methylprednisone, Hydralazine, Vancomycin, Ciprofloxacin, Fluconazole, D50, Insulin, Levothyroxine, Mannitol, Diuril, Nitroglycerin, Dopamine, Normal Saline, 20 KCL, Nipride, Vasopressin and Heparin. Flushed and stored in HTK
HMC309	Male	25	C	Head Trauma-MVA	Chewing Tobacco- Once monthly X 5 yrs. Beer- 4 per night daily. Liquor 2-3 drinks per night. Either or but not both in a given night.		Insulin, IVF, Ancef, Fentanyl, NS, Albuterol, Cozaar and Heparin. Flushed and stored in UW solution.
*HMC310	Male	41	C	Head Trauma-Blunt Injury	Cigarettes- 1/2 PPD since age 15. Beer- 1-2/Weekdays, 9-10/Weekends X 10 years.	COPD- Diagnosed within last 2 yrs. Pneumonia and Bronchitis within last yr. Hypertension; Diagnosed 2 yrs ago. Atenolol, Zyrtec, and Albuterol Inhaler (as per NOK).	Nicardipine, Insulin, Crystalloids, Albuterol, Clindamycin, Gentamycin, Keppra, Rocephin, and Zosyn. Flushed and Stored in HTK.
HFC312	Female	45	C	Anoxia	Cigarettes- 2.5 PPD X 22 yrs. Vodka- Heavy ETOH use. 1 Quart/ 3 days. Cocaine- Snorted- Used within the last 6 months.	Asthma/COPD- Pt used inhaler. Theophylline, Nasonex, Advair, Cymbalta, Neurotin, Albuterol, Chantix, and Nicoderm Patch(as per NOK)	Neosynephrine, Solumedrol, Insulin, IV Dextrose, Rocephin, Zithromax, Amiodorone, Heparin and Zemuron. Amiodorone, 4 rounds of Epinephrine, 2 rounds of Atropine, Narcan, Magnesium Sulfate administered during first response.
*HFC317	Female	46	C	Anoxia Anoxia-2nd to Cardiopulmonary Arrest	Cigarettes: 1/2-1 PPD X 26-30 yrs. Wine Cooler: 1 Every 2 weeks or less since age. Brandy/beer 4-6 times a yr X 25-30 yrs.	Emphysema- Diagnosed within 2 yrs of death. COPD- Diagnosed within 2 yrs of death. Hypertension- Tx w/Meds. Type II Diabetes: Diagnosed > 15 yrs prior to death. Insulin dependent X 10-15 yrs. Pt told could have fatty liver disease 5 yrs prior to death. Pt had bladder infections in year prior to death. Depression Meds, Heart Meds, Diabetes Meds, Insulin and Anxiety Meds (as per NOK). Admitted for CHF with weight gain of 20 pounds one week prior.	Neosynephrine, Epinephrine, Levophed, Vasopressin, Platelets, Regitine and Heparin. Flushed and stored in UW solution
*HFC336	Female	45	C	CVA	Beer- weekly	No Liver disease but in 1996 pt became jaundiced and a growth was found under the liver. It was removed, tested and determined not to be cancerous. No follow up required and no problems since the surgery. No home meds reported by NOK.	annitol, Hydralazine, Insulin, Heparin Sodium, NS Bolus, D50, T4, Zosyn, K-Phos, Lasix, Insulin, Methylprednisone, Fentanyl, Propofol, Dopamine, Levophed, Nicardipine, Vasopressin and Solumedrol. Flushed and stored in HTK.

For Research Use Only. Not for use in diagnostic or therapeutic procedures.

© Discovery Life Sciences

Discovery Life Sciences
 6 Henshaw Street
 Woburn, MA 01801
 Tel: (866) 838-2798
info@dls.com
<https://www.dls.com/>

UltraPool™ HLM 150

Donor Information

Specimen	Gender	Age (years)	Race	Cause of Death	Social History	Medical History	Medication given during Hospitalization
*HFC343	Female	65	C	ICH-Stroke	Cigarettes: 1 PPD X 17 yrs. Quit in 1979.	Asthma- tx with Albuterol. Grand Mal Seizures- Diagnosed in 1980. Recently diagnosed with Diabetes. Dilantin, Darvacet, Synthroid, ASA, Lasix, Albuterol, Enalapril, Pravastatin, Naproxen, and Hoodia (as per NOK)	Atropine, Dopamine, K rider, Lidocaine, Lasix, Solumedrol, K-Phos, Magnesium, Ancef, DDAVP, Insulin, Heparin and Nimbox. Epinephrine, Atropine, Lidocaine given at time of CPR. Flushed and stored in UW.
*HFC345	Female	69	C	Anoxia-2 nd to Cardiac Arrest	Wine- Socially.	Hypertension- diagnosed 4 months prior. Compliant with meds. Vitamin and HTN meds. (as per NOK)	Levophed, Dilantin, Insulin, KCL, T4, Vasopressin, Dopamine, Solu-Medrol, Calcium Gluconate, D5W, Albumin, Lasix, Heparin, Mannitol and NS. Flushed and stored in UW solution.
HMC349	Male	75	C	Hemorrhagic Bleed	Cigarettes; Quit in early 1990	Angina, CHF. Hypertension; Unknown length of time. Tx with Digoxin. Kidney Stones; early 1960's.	Vitamin K, Fee H2O, Cardizem, Zosyn, Mannitol, Nexium, 1/2 NS, Levophed, Vasopressin, Nipride, Heparin, and Vasopressors. Flushed and stored in UW solution.
HFC350	Female	54	C	ICH-Stroke	Not available.	Multi Vitamins and Calcium (as per NOK). No significant PMH.	Levophed, Levothyroxine, 1/4 NS, Neosynephrine, Lasix, Magnesium Sulfate, Calcium Chloride, Calcium Gluconate, NaHCO3, KPhos, KCL, Vitamin K, Ancef, Solumedrol, Insulin, Heparin, Vasopressors and D50. Flushed and stored in UW solution
HFC359	Female	60	C	Acute Respiratory Failure	Non-smoker; Non-drinker	Asthma- Not known when diagnosed. Hypertension- diagnosed 5 yrs prior. Controlled w/ Diovan. Diabetes- Diagnosed 6 yrs prior. Treated with Metformin. Hyperlipidemia- Diagnosed < 5 yrs prior. Controlled with Pravachol. GERD- Diagnosed 5 yrs prior. Treated with Prilosec. Prozac, Diovan, Klonopin, Prilosec, Pravachol, Metformin, Aspirin, Nasonex, and lycra (as per NOK)	Levophed, Solumedrol, Insulin, Ativan, Fentanyl and Morphine. Flushed and stored in HTK.
*HFC365	Female	42	C	Anoxia- 2 nd to suicidal drug overdose	Cigarettes- 1.5 PPD X 15 yrs. Quit 6 months ago. Vodka- 1 pint/week X 7 yrs. Quit in 2007. Benzodiazepine and opioid use.	Tizandine, Lortab, Rocaltrol, Trazadone, Ferrous Sulfate, Estradiol, Reglan, Januvia, Ambien, Phenergan, symbicort, Lamictal, Valium, Prozac, Remeron, and Lortab (as per NOK)	Dopamine, Keppra, Doripenem, Linezolid, Insulin, Clindamycin and Heparin. Flushed and stored in UW Solution.
HFC372	Female	73	C	ICH-Stroke	Cigarettes- 2 PPMonth X 17 yrs. Quit 20 yrs. Ago.	Cardiac Disease- NOK unsure of diagnosis. Hypertension- Treated with meds. Diverticulitis. Treated with meds. Dicyclomine, Atenolol, Plavix, Diovan, Simvastin, Hyomax, Prednisone, Omeprazole, ASA, Synthroid, and Xanax (as per NOK)	Levophed, Neosynephrine, Vasopressin, Dopamine, Vancomycin, Crystalloids, Colloids, Heparin, Mannitol and Lasix. Flushed and stored in UW solution.

For Research Use Only. Not for use in diagnostic or therapeutic procedures.

© Discovery Life Sciences

Discovery Life Sciences
6 Henshaw Street
Woburn, MA 01801
Tel: (866) 838-2798
info@dls.com
<https://www.dls.com/>

UltraPool™ HLM 150

Donor Information

Specimen	Gender	Age (years)	Race	Cause of Death	Social History	Medical History	Medication given during Hospitalization
*HFC377	Female	54	C	CVA-ICH	Cigarettes- 1 PPD X 30 yrs. Quit in Jan 2005.	Bypass surgery- 1999. Hypertension- Diagnosed 8 yrs prior. Nitro (as per NOK)	Neosynephrine, Zosyn, Solumedrol, Albuterol, T4, Insulin, Dopamine, Adenosine, Bicarb, 5% Albumin, ICA, Levophed, FFP, PRBCs, Crystalloids and Heparin (administered in hospital) Flushing Solution: UW. Storage Solution: UW.
*HFC379	Female	37	C	ICH-Stroke	Cigarettes; 2-3 PPD x 10-15 yrs. Vodka/wine coolers- on occasion. Marijuana- experimented with in HS.	Hypertension; diagnosed in early 20's. Xanax, BP Meds, and Multi Vitamin (as per NOK).	Dopamine, Levophed, Amiodarone, Heparin, Mannitol, Vasopressors and Crystalloids. Flushed and stored in UW solution.
*HFC380	Female	45	C	Anoxia-2 nd to Respiratory Failure	Non-smoker; Non-drinker	Severe Asthma >30 yrs. Has been hospitalized over 10 times for Asthma related reasons.	Levophed, Fentanyl, Cardene, Ativan, Crystalloids, Vasopressin, Zemuron, Combivent, Fragmin, Insulin, Lasix, Levothyroxine, Mannitol, Narcan, Solumedrol, Zosyn, Zyvox, NS, Pitressin, T4, Heparin and Vasopressin. Flushed and stored in HTK.
*HFC430	Female	50	C	Anoxia-2 nd to Cardiovascular	Cigarettes- 2 PPD X 25 yrs. Vodka & Mixed Drinks- Daily X 25 yrs. Cocaine & Marijuana- 6 yrs ago.	Asthma- Developed as an adult. Hypertension- Diagnosed 10 yrs prior. Elevated liver enzymes upon admit. HCTZ (as per NOK)	Propofol, Unasyn, D5 1/2 NS, KCL, Metoprolol, Heparin and Thorazine. Flushed and stored in HTK.
*HFC449	Female	57	C	ICH-Stroke	Non-smoker; Non-drinker	Hypertension- Diagnosed 3 yrs prior. Compliant with meds. Recent Seizure upon admit. Type II Diabetes- Diagnosed 6 yrs prior. Compliant with meds. Zydis, Lovastatin, Acetaminophen, Aspirin, Bupropion, Dicyclomine, Escitalopram, Flurazepam, Gabapentin, Lisinopril, Metformin, Ziprasidone (as per NOK).	NS, Dopamine, Levaquin, Clindamycin, Insulin, Levothyroxine, Narcan, Norcuron, Pitressin, Potassium Chloride, Sodium Bicarbonate, Solumedrol, Water, Albuterol, Atrovent, Magnesium.
HFC452	Female	64	C	ICH-Stroke	Wine- 1 glass every 3 days X 3-4 yrs.	Urinary Tract infection- 2009. Belbutrain, Calcium, and Magnesium Phosphate (as per NOK).	Levophed, I.V. Fluids, Solumedrol, Lasix, Arginine Vasopressin, PRBCs, Hespan and T4. Flushed and stored in UW solution.

For Research Use Only. Not for use in diagnostic or therapeutic procedures.

© Discovery Life Sciences

Discovery Life Sciences
 6 Henshaw Street
 Woburn, MA 01801
 Tel: (866) 838-2798
info@dls.com
<https://www.dls.com/>

UltraPool™ HLM 150

Donor Information

Specimen	Gender	Age (years)	Race	Cause of Death	Social History	Medical History	Medication given during Hospitalization
*HFC458	Female	44	C	Anoxia	Cigarettes- 2 PPD X 30 yrs. Patient smoked marijuana as teenager- Not for long.	Hypertension. Metoprolol, Lisinopril, Prozac, Ambien, Percocet, and Xanax (as per NOK). Significant depression and heart palpitations, secondary to anxiety	Sodium Bicarbonate, Zosyn, Potassium Chloride (k+), Heparin, Potassium Phosphate (K-Phos), Calcium gluconate, DDAVP, D50, Methylprednisone, NS, Naloxone, Levothyroxine, Vasopressin, Albuterol, Levophed, Insulin, Dopamine, 0.45 NS, Levothyroxine, D5W, Vasopressin, PRBC, 5% Albumin, KCl in NS, D5, Mannitol, T4, and Crystalloid. Flushed and stored in UW solution
*HFC465	Female	66	C	ICH-Stroke	Cigarettes- Smoked for 10 yrs. Quit in 1972. Wine with dinner. Occasional ETOH.	Asthma with chronic cough. Carried emergency inhaler. Baby Aspirin, Estrogen Patches, Nortriptyline, and Prilosec (as per NOK).	Levophed, Dopamine, Amiodarone, Vassopressin, PRBCs, Platelets, D5 1/4 NS, FFP, Insulin, Albumin 5%, Lasix, Solumedrol, Ampicillin-Sulbactam, Dexamethasone, Crystalloids, and Heparin. Flushed and stored in UW solution.
HFC469	Female	62	C	ICH-Stroke	Cigarettes: 1 PPD X 10 yrs. Wine: Occasional glass X 40 yrs.	COPD: Diagnosed 5 yrs prior. A-Fib, pacemaker: 2 yrs. Diagnosed with Hypertension 15 yrs prior. Basal cell on nose: Removed 4 yrs ago. Diagnosed with Diabetes 15 yrs prior. Insulin dependent for 10 years. Insulin and HTN Meds (as per NOK)	Norepinephrine, Levothyroxine, T4, Magnesium Sulfate, Digoxin, Docusate, Furosemide, Colace, Isosorbide, Omeprazole, Tylenol, Hydralazine, Labetolol, Cefapriime, Vancomycin, Potassium, Phosphate, Glycolpyrrolate, Insulin, Hetastarch, Heparin, and Levophed. Flushed an stored in HTK.
HFC474	Female	52	C	ICH-Stroke	Cigarettes: 1.5 PPD X 28 yrs. Quit 4 yrs ago. Drank Socially.	Diagnosed with Hypertension 1 yr prior. Unknown HTN Med (as per NOK)	Amiodarone, Phenylephrine, Mannitol, Metoprolol, Albuterol, NS, Zosyn, Neosynephrine, Propofol, NaPhos, Vancomycin, Crystalloids, Colloids, Heparin and Solumedrol. Flushed and stored in UW.
*HMC487	Male	54	C	Anoxia-2 nd to Cardiovascular	Cigarettes: 1 PPW X 20 yrs. Quit many yrs ago. Wine: as a young man. Marijuana: Smoked yrs ago. Cocaine: Snorted 15 yrs ago. Shrooms: Used yrs ago.	Dementia 2 nd to Parkinson's Disease within last 10 yrs. Many (as per NOK).	Xopenex, Ceftriaxone, Lovenox, Famotidine, Methylprednisone, Cefazolin, 1/2 NS, Kphos, Ancef, and Heparin. Flushed and stored in UW.

For Research Use Only. Not for use in diagnostic or therapeutic procedures.

© Discovery Life Sciences

Discovery Life Sciences
 6 Henshaw Street
 Woburn, MA 01801
 Tel: (866) 838-2798
info@dls.com
<https://www.dls.com/>

UltraPool™ HLM 150

Donor Information

Specimen	Gender	Age (years)	Race	Cause of Death	Social History	Medical History	Medication given during Hospitalization
*HFC488	Female	43	C	Anoxia-2 nd to Cardiac Arrest	Cigarettes: 4-5 ciggs X 3 yrs. Beer: 1-2 rarely X 25 yrs.	Normal healthy	Fentanyl, Sodium BiPhos, Glutamine, Protonix, Lovenox, Oxycodone, Ampicillin, Chloraseptic, Vancomycin, Zosyn, Tylenol, Motrin, Sodium Chloride, Insulin, RBCs, FFP, Dopamine and Levophed. Flushed and stored in LR/UW
HMC489	Male	82	C	CVA-Stroke	Non-smoker; Non-drinker	Hypertension: Diagnosed 10-15 yrs ago. Type II Diabetes: Diagnosed last October. Bleeding Ulcers: 1953- No reoccurrence. Avodor, Glemepeside, Simvastin, Enapril/HCTZ and Lubiprostone (as per NOK).	T4, Neosynephrine, Crystalloids, Ancef, Heparin, Mannitol, and Lasix. Flushed and stored in UW solution.
*HMC490	Male	77	C	CVA-ICH	Non-smoker; Non-drinker	Bronchiectasis- lung labectomy in 1956. Used SVN machine twice a day. 2 stents placed. Hypertension: Diagnosed 25 yrs ago. Prostate Cancer: Diagnosed 13+ yrs ago. Surgically removed and resolved with no other treatment. Colon surgery 3 yrs ago to remove dead tissue caused by HTN. Aspirin, Plavix, Lipitor, Diovan, and Amlodipine (as per NOK).	Solumedrol, T4, Insulin, Arginine Vasopressin, and Vecuroinum. Flushed and stored in HTK.
*HFC492	Female	41	C	ICH-Stroke	Wine: Occasional glass.	CHF: Diagnosed 2 yrs prior. Stent placed. 3 Heart Caths and an MI. History of CADD and A-fib. Hypertension: Diagnosed 15 yrs prior. Type I Diabetes since age 11. Treated with Insulin injections. GERD & Gastritis. Hernia repair surgery. Chronic Kidney Disease. Kidney/Pancreas tx in 1195- rejected both 11 days later. 2nd Kidney/Pancreas tx in 2000- rejected pancreas.	Acyclovir, Amlodipine, Benzonatate, Ciprofloxacin, Clopidogrel, Cyclosporine, Aranesp, Albumin Free, Ferrous Sulfate, Folic Acid, Furosemide, Glucagon, rDNA, Hydralazine, Novalog, Humulin, Imdur, Labetolol, Cellcept, Nitroglycerin, Omeprazole, Prednisone and Simvastin. Labetolol, Dopamine, Norepinephrine, Levothyroxine, Vasopressin, Insulin, Normal Saline, Potassium Chloride, Albumin, Sodium Bicarbonate, Lasix, Dextrose, Methylprednisone and Heparin. Flushed and stored in HTK
*HFC494	Female	55	C	CVA-Stroke	Cigarettes: 1.5 PPD X 8 yrs. Marijuana: Occasional smoker.	Hypertensive from 1998-2003- treated with meds. Resolved after gastric bypass surgery in 2003. Seizures: Diagnosed 4-5 yrs prior. 1-3 times a month as of lately. Diverticulitis: Diagnosed before 2003. Dilantin, Keppra, Paxil, Nexium, Fentanyl and Oxycodone (as reported by NOK)	Insulin, PRBCs, Heparin, and Solumedrol. Atropine and Epinephrine administered at time of CPR. Flushed and stored in UW.

For Research Use Only. Not for use in diagnostic or therapeutic procedures.

© Discovery Life Sciences

Discovery Life Sciences
6 Henshaw Street
Woburn, MA 01801
Tel: (866) 838-2798
info@dls.com
<https://www.dls.com/>

UltraPool™ HLM 150

Donor Information

Specimen	Gender	Age (years)	Race	Cause of Death	Social History	Medical History	Medication given during Hospitalization
*HFC499	Female	28	C	Head Trauma-2 nd to MVA	Cigarettes: 2.5 PPD X 13 yrs. Beer: Bottle a day on weekends. Marijuana: Smoked every 3 months X 10 yrs. Last used a week prior.	Mild Bronchitis- 2-3 times a year. Treated with Duoneb. Pre Clampsia- With both pregnancies. Last time 6 months. Fish Oil (as reported by NOK)	V Dextrose, Insulin, PRBCs,
*HMC500	Male	54	C	CVA-ICH	Beer: Occasionally 1-2 times a week since 18 yrs old.	CVA- 3 months ago requiring 1 wk hospital stay. CVA: 7 yrs ago. Lexapro, Lipitor, ASA, Altace, Coumadin, and Alovert (as per NOK).	Crystalloids, Colloids, Cardene, Vancomycin, Platelets, and Heparin. Flushed and stored in UW solution.
HMC501	Male	76	C	ICH-Stroke	Cigarettes: 2 PPD X 30 yrs. Quit in 1993. Whiskey: 8 shots per wk x 15 yrs. Quit November 2010.	Congestive Heart Failure- Treated with BP meds and water pills. Hypertension: Diagnosed 30 yrs ago. Compliant with meds. Multivitamin and aspirin (as per NOK).	Labetalol, NS, FFP/ Plasma, Vitamin K, Insulin, Levothyroxine, Potassium Phosphate, Solumedrol, KCL, Zosyn, Nicardipine, Heparin, and Zemuron. Flushed and stored in HTK.
*HFC506	Female	72	C	Anoxia-2 nd to Cardiovascular	Social alcohol at holidays X adult life.	History of Asthma- Treated with Advair. Hypertension- Diagnosed > 40 yrs ago. Treated with Norvasc. History of Gait changes- Vertigo. Type II Diabetes X 15 yrs. Insulin dependent X 15 yrs. Irritable bowel syndrome. Kidney Failure X 4 yrs. Kidney Stones over last couple years. Treated with Kidney Dialysis in 2001. Frequent bladder infections & UTIs. Norvasc, Advair, Acetaminophen, Aspirin, Florajen, Gemfibrozil, Humalog, Lantus, Percocet, Prednisone, Prograf, Bactrim and Toprol as per NOK).	IV Fluids, Methylprednisone, Hydrocortisone, Insulin, Vasopressin, Dopamine, Levophed, Bicarb, Heparin, Calcium, Invanz, Lasix, Albumin, Magnesium, Kphos, KCL, Azithromycin, Vancomycin, Meropenum, Pitressin, CaCl, LR, 0.9 NS, and Hespán. Flushed and stored in UW solution.
*HFC507	Female	61	C	ICH-Stroke	Cigarettes- Smoked < 1 Pack per yr socially. Alcohol use very infrequently: once a year.	History of CHF, Pacemaker, and A-Fib. Hypertension- Diagnosed 35 years- Compliant with meds. Diabetes- Type II X 2-3 yrs. Took oral meds & controlled with diet. Gabapentin, Lisinopril, Januvia, Metoprolol, Metolazone, Hydralazine, Bumetanide, Klor Con, Isosorbide, Coumadin, ASA, and Vege Lax (as per NOK).	Dopamine, IV Fluids, Lasix, Bumex, T4, FFP, Bicarb, Neosynephrine, Levophed, Dobutamine, Epinephrine, and Heparin. Flushed and stored in UW solution.
*HMC509	Male	50	C	Complications-2 nd to Lupus	Beer; 1-2 per day socially x past few years & short time when he was young. Marijuana: Smoked frequently/few times per day x 15 yrs and short time when he was young.	Nodules found in lungs. Rare use of inhaler. Calcified Valves. Aortic stenosis; recently diagnosed. Hypertension; diagnosed 10-12 yrs ago. Esophageal Ulcer; treated with meds & diet. Kidney Stones; 16 yrs prior & 7 yrs prior. Meds for stomach, Heart Meds, ASA, Lopressor, HTN meds, Altace, and Valium (as per NOK).	IV Fluids, Dextrose, Milranone, Levophed, Hydrocortisone, Prednisone, Lasix, Insulin, Vasodilators, Arginine Vasopressin, PRBCs, FFP, Cryo, Platelets, Albumin.
*HMC510	Male	22	C	Head Trauma-2 nd to Blunt Injury	Beer and Liquor: Socially since 18 yrs old.	Healthy	Levophed, IV Fluids, Insulin, Arginine Vasopressin, Pentobarbital, Fentanyl, Heparin, Dobutamine, and Ancef. Flushed and stored in UW solution.

For Research Use Only. Not for use in diagnostic or therapeutic procedures.

© Discovery Life Sciences

Discovery Life Sciences
 6 Henshaw Street
 Woburn, MA 01801
 Tel: (866) 838-2798
info@dls.com
<https://www.dls.com/>

UltraPool™ HLM 150

Donor Information

Specimen	Gender	Age (years)	Race	Cause of Death	Social History	Medical History	Medication given during Hospitalization
*HMC513	Male	68	C	ICH-Stroke	Cigarettes: 1 PPD since age 17. Quit in 1986. Beer: 1 a week socially since a teenager.	COPD: Treated with meds due to heart problems. CAD/CHF since 1987. CABG X 2. Stents. History of LAVD. Hypertension: Diagnosed in 1987. Diabetes in 1987. Treated with Novolog and Lantis. Insulin dependent X 20 yrs. Meds for HTN, Diabetes, Depression, Heart, Multivitamin, Magnesium, Folic Acid and Protonics (as per NOK). History of LVAD and Von Willebrand Disease.	Propofol, Heparin, Coumadin, Insulin, Levophed, amitriptyline, Aspirin, Carvediol, Cetrizine, Divalproex, Docusate, Folic Acid, Hydrazaline, Hydrocodone, Insulin Glargine, Insulin Lispro, Isosorbide, Lisinopril, Magnesium Oxide, Magnesium Sulphate, 25% Mannitol, Methimazole, Morphine, Multivitamin, Naloxone, Pantoprazole, Potassium Chloride, Prochlorperazine, Rosuvastatin, Topiramate, Venlafaxine, Warfarin Sodium, Spironolactone, FFP, 5% Albumin, NS, Pitressin, Narcan, Solumedrol, KCL, Kphos, Ancef, Heparin, and Hespan. Flushed and stored in UW solution.
*HFC518	Female	42	C	CVA 2 nd ICH	Non-smoker; Non-drinker	Donor was prescribed Paxil; reason not mentioned in NOK questionnaire. Donor took antidepressant Paxil, Ambien as needed and not sure if still taking cholesterol medication (as per NOK).	Neosynephrine, Levophed, Insulin, IVPB, Vasopressin, 1/2 NS, D5 1/4 NS/20kcl, 1/2 NS/bicarb, KCL, Lasix, DDAVP, Albuterol, Unasyn, Levothyroxine, Cefalozin, Solumedrol, Levaquin, Narcan, Heparin, Crystalloids, Epinephrine, Norepinephrine, and Dopamine, Flushed and stored in SPS.
*HFC519	Female	54	C	CVA 2 nd ICH	Non-smoker; Non-drinker	Donor had no known cancer, however, suffered with benign ovarian cysts which led to a hysterectomy. Donors chart shows "insulin resistance" and DM type 2. NOK states none in medical/social history. Donors medical/social history states NO liver disease, however, interviewer mentions in notes further complications of liver biopsy in 2009. Donor was diagnosed with Celiac Disease in 2009. Lisinopril, Zanaflex, Remeron, Urodiol, Ritalin, Klor-con, Klonopin, Levothyroxine, Lasix, Trazadone, and Nitrofurantoin (as prescribed by physician).	Levophed, Bicarb, Xigris, Solumedrol, insulin, Vecuronium, Zosyn, Zithromax Succinylchloride, Etomidate, Narcan, Rocephin, Heparin and Crystalloids. Flushed and stored in UW solution.
HMC521	Male	25	C	SIGSW	Abuse rehab 7 years ago; alcohol, cocaine (inhaled). Positive for alcohol on admission.	Herpes.	Dopamine, D5, KCl, potassium, Ancef, DDAVP, Hespan.

For Research Use Only. Not for use in diagnostic or therapeutic procedures.

© Discovery Life Sciences

Discovery Life Sciences
6 Henshaw Street
Woburn, MA 01801
Tel: (866) 838-2798
info@dls.com
<https://www.dls.com/>

UltraPool™ HLM 150

Donor Information

Specimen	Gender	Age (years)	Race	Cause of Death	Social History	Medical History	Medication given during Hospitalization
*HFC523	Female	51	C	Anoxia-2 nd to Drug Intoxication	Cigarettes- 1/2 PPD X 30 yrs. Beer- 6-8 beers on weekends. Marijuana- Smoked once in a while over past 30 yrs.	Diagnosed with respiratory disease several yrs ago. Treated with Albuterol. Diagnosed Hypertensive 2 years prior. Med compliant. Norvasc, Albuterol, Lopressor, Fexofenadine, Multivitamin, ASA, Wellbutrin, Trileptal, Fish oil, Vitamin C and Lisinopril (as per NOK). Pt is believed to have taken Restoril, Norvasc, Klonopin and Elavil.	Norepinephrine, Dopamine, Vasopressin, Zosyn, KCL, Calcium Gluconate, Solu-Cortef, DDAVP, Insulin, D5W, and Heparin. Flushed and stored in HTK.
HMC526	Male	48	C	Anoxia	No tobacco or drug use.	Diabetes, seizures, obesity.	Labetalol, Nipride, Lidocaine.
HFC529	Female	64	C	CVA 2 nd ICH	Pt smoked for 3-4 yrs during college but quit after graduating and has not smoked since. Alcohol; Seldom and on special occasions.	Pituitary tumor; Diagnosed 15-20 yrs ago. It was found to be benign and checked annually. Pt took Vitamins and a calcium supplement daily (as per NOK).	Pitressin, D5, Nicardipine, Insulin, Narcan, Hydralazine, NS, Ancef, Norcuron, Atrovent, Albuterol, Humulin, Cardizem, Vecuronium, Neosynephrine, Solumedrol, Insulin, Vasopressin, Heparin, Antihypertensives. Flushed and stored in HTK.
HMC530	Male	30	C	Intracranial Bleed	Non-smoker, occasional drinker	No medical history available	Dopamine
HMC634	Male	57	C	Subarachnoid hemorrhage	Whiskey.	Heart disease >10 years. Arthritis, stiff joints. Hernia operation 13 years ago. Arthroscopic knee surgery 5 years ago. Indoral, HCTZ, aspirin. Amoxicillin x2 weeks.	Heparin, Epinephrine, Zemuron, Dopamine, Hespan, Piperacillin, Cipro, Diflucan, crystalloids, colloids.
*HMC657	Male	64	C	Intracranial Bleed	Heavy alcohol use in past, cut down in last year.	Hypertension. Thyroid problems. Asbestos exposure 30 years ago. Blood in urine 2 weeks ago; renal carcinoma noted on kidney biopsy. Synthroid, Glynase, Ceflex.	Nipride, Deiprivan, Decadron, Pepcid, Dilantin, Nitro paste, Nimodipine, Ancef, KCl, Mg
*HMC712	Male	45	C	Head Trauma	Smoker, Infrequent vodka drinker	Cluster migraines. Elevated blood pressure 2nd to anxiety. Substance abuse problem last 10 yrs for Tylenol w/codeine, hydrocodine, Percocet off/on. Daily meds Anacin, Aleve, Prozac, Catapress, Tylenol w/codeine, Percocet	Dopamine, Mannitol, Decdron, Zantac, Pitressen, Succinylcholine
*HFC726	Female	71	C	ICH	3 or 4 drinks a week.	History of arteriovenous malformations. Prior ICBs 1965, 1967, 1976. Knee replacement 1997. Mitral valve replacement 1995. Hemorrhoids. Skin cancer on nose 5 or 6 years ago; malignant, not melanoma. Cataract surgery 2-3 years ago. Coumadin, Detrol, Celebrex.	Heparin, Dopamine, KCl, Ancef, insulin, D5, Vit K, blood products.
*HFC738	Female	55	C	Cerebrovascular Accident, Intracranial hemorrhage	Normal.	DM x8 years, IDDM x2 years. Hypertension, CAD x20 years, angioplasty 1991. MI x2. Cardiac catheterization 1996. CVAs, TIAs x6 years. Endarterectomy 4 weeks prior. D&C x2, hysterectomy 10 years ago. Glaucoma x2 weeks. Appendectomy, tonsillectomy as child. Insulin, aspirin, Zocor, Glipizide, Lopressor, Isordi, Persantinel. Vaccinated for Hep B.	Heparin, Lasix, Dopamine, KCl, blood products, crystalloids.

For Research Use Only. Not for use in diagnostic or therapeutic procedures.

© Discovery Life Sciences

Discovery Life Sciences
6 Henshaw Street
Woburn, MA 01801
Tel: (866) 838-2798
info@dls.com
<https://www.dls.com/>

UltraPool™ HLM 150

Donor Information

Specimen	Gender	Age (years)	Race	Cause of Death	Social History	Medical History	Medication given during Hospitalization
*HMC778	Male	65	C	Cerebrovascular Accident	History of alcohol abuse; quit drinking 1972. Possible jail time for drunk and disorderly as youth.	Endocarditis; aortic valve replacement 1987. Rotator cuff surgery 1981, disc surgery 1982, eye surgery 1980. Kidney stones 1998. Hypertension x20 years, controlled by meds. Diabetes x1year, controlled with diet. Mild asthma, no medications. Coumadin, Cardizem, KCl, Anticard. Concussion week prior to admission - kicked in head by horse.	Norcuron, Lasix, Mannitol, Vit K, Dopamine, LCNW, Vasopressin, Insulin, T4, Nipride, blood products, crystalloids
*HMC786	Male	22	C	Head trauma, MVA	Social alcohol. Positive for marijuana on admission, no social history of drug use.	Biopsy on thyroid nodule 10/01, results negative.	Heparin, Dopamine, crystalloids, tetanus shot, Cefazolin, Clindamycin, Ancef, albumin, Zantac, Mylanta, DDAVP, KCl
HMC800	Male	33	C	Head trauma, ped v. MVA	6 pack beer per day, unknown duration. Marijuana, unknown duration. Multiple tattoos.	Broke right collar bone 2 years ago. Foreign body in right foot, 3 years, never removed. MVA 2 months ago. Allergic to topical ointments. Dog bite 3 years ago; no symptoms.	Heparin, blood products.
HFC839	Female	49	C	Anoxia s/p cardiovascular	None available (no next of kin)	Downs Syndrome. Valve disease, mitral stenosis, endocarditis secondary to AV graft infection 7/04 with MSSA and pseudomonas. Pulmonary hypertension, CRF with hemodialysis, diabetes. A Fib, hypertension, colitis. Coumadin. Grand mal seizure during hospitalization, nothing noted on neuro history.	Heparin, Mannitol, Dobutamine, Neosynephrine, Levophed, Dobutex, Dopamine, Dilantin, Cefepime, Solumedrol, Humulin, Topramycin, Flagyl. (records unclear)
*HFC851	Female	39	C	ICB s/p MVA	Drank liquor and beer daily, large quantities. Cocaine use, unknown route.	Caesarean section, tubal ligation 2004; varicose veins.	Heparin, Neosynephrine, insulin, Phenobarbital, Atracurium, Mannitol, Labetalol, Pepcid, Ancef, Dopamine, Remifentanyl
*HFC903	Female	60	C	Anoxia-2 nd to Seizure	Drank 1 glass wine socially X 15 years.	Hypertension- Diagnosed in 2010. Compliant with meds. Seizures from Epilepsy since age 7. Seizure meds, HTN meds, Plavix, and multivitamin (as per NOK). Plavix, Lamictal, Naproxene, Crestor, Topamax, Bystolec, Maxide, Lopressor, and Topyramaz Lvyxyl.(taken at home per OPO donor form).	Zosyn, Solumedrol, Albuterol, Primaxin, Keppra, Lopressor, Cefapime, 1/2 NS, Heparin, PRBC 250 and FFP 250. Flushed and stored in UW solution.
*HFC908	Female	70	C	Anoxia 2 nd to Cardiovascular	Wine; Daily.	COPD/Pneumonia July previous yr and currently. Colon polyps removed 10 yr prior. Actonel, Armour Thyroid, Calcium Folic Acid, Methotrexate, Mobic, Multivitamin with minerals and Remicade (as per NOK).	Dopamine, NS, KCL, Levothyroxine, Bicarb, Vasopressin, Ancef, Vitamin K, Solumedrol, Insulin, D50, Armour Thyroid, Lidocaine, Albuterol, Elavil, Fentanyl, NaCl, Crystalloids, T4, Vecuronium, Heparin, and Lasix. Flushed and stored in UW solution.

For Research Use Only. Not for use in diagnostic or therapeutic procedures.

© Discovery Life Sciences

Discovery Life Sciences
 6 Henshaw Street
 Woburn, MA 01801
 Tel: (866) 838-2798
info@dls.com
<https://www.dls.com/>

UltraPool™ HLM 150

Donor Information

Specimen	Gender	Age (years)	Race	Cause of Death	Social History	Medical History	Medication given during Hospitalization
*HFC909	Female	64	C	Anoxia 2 nd to Cardiovascular	Cigarettes: smoke 2 PPD since 15 yrs old. Beer: Drank unknown amount since 20 yrs old, frequency unknown (per NOK).	COPD; Donor suffered with COPD since 1 yr ago. Hypertension: Donor suffered for 4 yrs and it was controlled by diet and exercise. Donor did not take any prescribed medications (as per NOK).	Epinephrine, Atropine, Levophed, Amiodarone, Solumedrol, Ancef, Potassium, Albumin, Pepcid, Lovenox, Zofran, Mannitol, Morphine, Insulin and Aspirin. Flushed and stored in UW solution.
*HMC911	Male	53	C	CVA 2 nd ICH	Cigarettes; 1 PPD x 30-35 yrs; quit for 2 months 5 yrs ago. Whiskey and beer; pt drank occasionally for years; very little in the last 5 yrs. Marijuana; once per day daily x 38 yrs. Pt sniffed 3-4 lines of cocaine on special occasions, approximately once per year last used December 5 years prior.	Pt suffered Heart Valve prolapse 10 yrs prior. CVA. Hypertension; Pt suffered for 6 months; medication prescribed to control, but pt was non-compliant. Vertigo 5 yrs ago. Donor took Tums and Tylenol as needed and Lopressor for hypertension (as per NOK).	Neosynephrine, Dopamine, Mannitol, Solumedrol, Insulin, Heparin, Lasix, Thorazine, Crystalloids, T4, Nitroglycerine, and Vasopressin. Flushed and stored in UW solution.
HMC912	Male	41	C	Head Trauma-2 nd to MVA	Cigarettes; 3-4 cigs a day since age of 24. Beer; 2 beers every other day since age 14. Tequila; 2 shots every other day since age 14. Marijuana; Smoked since age 14.	Asthma as a child; outgrown as an adult.	Versed, MGSO ₄ , and Heparin. Flushed and stored in UW solution.
*HMC913	Male	37	C	Anoxia 2 nd to Asphyxiation	Beer: 2 per week.	Respiratory disease: Treated for TB with pills X 6 months in 1995. Melanoma- removed small sample within last yr- recommended to remove larger sample- pt never followed up. Jaundice at birth. Suspected IBS since 2000. Anti-depressants (non compliant per NOK).	Insulin, Levophed, Labetolol, Lasix, Zosyn, Vancomycin, Levaquin, Crystalloids, Colloids, NaCL, D5NS, Heparin, and Solumedrol. Flushed and stored in UW solution.
HFC915	Female	59	C	Head Trauma- Blunt Injury	Non-smoker; Non-drinker	Vitamins, Loratidine as needed, and Tobramycin eye drops (as per NOK).	Neosynephrine, Levophed, Dopamine, IV fluids with Solucortef, T4, DDAVP, Arginine vasopressin, Heparin, Mannitol, and Zemuren. Flushed and stored in UW w Insulin.
HFC921	Female	33	C	Anoxia / Cardiovascular	Cigarettes, 1/2 PPD since age 15. Beer, "a couple socially per week" X 10 yrs but none in last yr.	Interstitial pneumonitis. Fentanyl, Oxycodone, "many others unknown names" (as per NOK).	IV Fluids with Solumedrol, Combivent, Prednisone, T4 and Insulin, Heparin, Mannitol, Lasix, and Ancef. Flushed and stored in UW solution.

For Research Use Only. Not for use in diagnostic or therapeutic procedures.

© Discovery Life Sciences

Discovery Life Sciences
6 Henshaw Street
Woburn, MA 01801
Tel: (866) 838-2798
info@dls.com
<https://www.dls.com/>

UltraPool™ HLM 150

Donor Information

Specimen	Gender	Age (years)	Race	Cause of Death	Social History	Medical History	Medication given during Hospitalization
HMC922	Male	42	C	Other 2 nd Seizures	Non-smoker; Non-drinker	Closed head Injury at age 9 and Craniectomy, HTN X 5 yrs; compliant w/ meds. Lennox-Gastuat (Seizure Disorder)and intractable Seizure Disorder; Unknown when diagnosed; Unknown if compliant w/ meds. Depakote, Ranitidine, Lactulose, Levothyroxine, Phenobarbital, Alendronate, Miralax, clobazam, Keppra (as reported in hospital chart).	Neosynephrine, Lasix, Vancomycin, Zosyn, Gentamicin, Cefepime, Pentobarbital, Lamotrigine, Enoxaperine, Phenobarbital, Lamictal, Floxacin, Miralax, Synthroid, Lovenox, Ativan. Flushed in Custodial. Stored in HTK.
*HFC924	Female	26	C	Head Trauma 2 nd Blunt Injury	Cigarettes- 1 PPD X 12 yrs. ETOH (Beer, Wine, Liquor)- < 1 drink/ mos. Methamphetamine- Unknown amount X 4 mos.; Unknown frequency.	Post-Partum Depression. Pristiq, Klonopin (as per NOK).	Lactated Ringer's, Magnesium Sulfate, Calcium Chloride, Solumedrol, Zosyn, Albuterol, Vitamin K, Levophed, Vasopressin. Flushed and stored in HTK.
*HFC925	Female	46	C	CVA 2 nd ICH	Cigarettes- 1 PPD X 11 yrs. ETOH (Beer)- 1 drink every 3 days X 11 yrs.	HTN x Unknown amount of years but its stated as several; compliant with meds. Antidepressants, Beta Blockers for HTN (as per NOK).	Levophed, Vasopressin, Ertapenem, Norepinephrine, Calcium Chloride, Magnesium Sulfate, Potassium Chloride, Calcium Gluconate, Sodium Phosphate, Zosyn, Heparin. Flushed and stored in HTK
HMC928	Male	37	C	ANOXIA 2 nd SI ASPHYXIATION	Cigarettes: 1 PPW X 19 yrs. ETOH (Beer): Binge drinker, would stop for months then drink 18 at once. Marijuana (Smoked) weekly from age 15 to 17. (as reported by NOK).	Healthy	Epinephrine, Lidocaine, Dextrose, Ciprofloxacin, Heparin, Normal Saline. Flushed and stored in Custodial.
HMC929	Male	57	C	CVA 2 nd ICH	Unknown if Cigarettes or other Tobacco products: 1 1/2 PPD X 30 yrs. Marijuana (Smoked); Unknown quantity or frequency; Crack Cocaine (Smoked); Unknown quantity or frequency; Hydrocodone; Unknown quantity or frequency; quit using all drugs 5 yrs ago.	CAP and Left sided Pneumonia. Legionella found in Urine. Retrocardiac left lower lobe infiltrate. Effusion not excluded. Stable position of line and tubes. 3. Pulmonary venous congestion. GERD; Unknown If compliant w/ meds. Simvastatin, Enalapril, Prilosec, Advil (as per NOK).	Dextrose, Solumedrol, Insulin, Heparin. Flushed and stored in SPS-1.
HMC930	Male	56	C	Anoxia 2 nd Electrocutted	ETOH (Beer) 6 Beers every weekend; Unknown duration.	HTN x 4 yrs; Compliant w/ meds. Kidney stones. Lopressor (as per NOK).	Heparin. Flushed and stored in SPS-1

For Research Use Only. Not for use in diagnostic or therapeutic procedures.

© Discovery Life Sciences

Discovery Life Sciences
 6 Henshaw Street
 Woburn, MA 01801
 Tel: (866) 838-2798
info@dls.com
<https://www.dls.com/>

UltraPool™ HLM 150

Donor Information

Specimen	Gender	Age (years)	Race	Cause of Death	Social History	Medical History	Medication given during Hospitalization
*HFC931	Female	19	C	Anoxia 2 nd Seizures	Non-smoker; Non-drinker	Seizures; Part of the current unexplained illness. No drug use (as per NOK).	Norepinephrine, Solumedrol, Insulin, Vasopressin, Heparin, Fentanyl, Pentobarbital, Lovenox, Magnesium Sulfate, Doxycycline, Propofol, Phenytoin, Midazolam, Ceftriaxone, Vancomycin, Acyclovir, Ancef, Calcium Gluconate, Levetiracetam, Potassium Chloride. Flushed and stored in HTK.
*HMC933	Male	39	C	Head Trauma 2 nd Blunt Injury	10 Cigarettes / weekend X 20 yrs. ETOH (Beer) 12-24 cans during the weekend X 20 yrs. Marijuana (smoked) Unknown amount used or frequency X 20 yrs; Pt had not smoked in last 3 months.	HTN; Diagnosed 5 yrs ago; unknown if compliant w/ meds. Kidney stones; Passed Kidney stone in 3 separate occasions. Vasodilators, Diuretics, Seasonal allergy meds, anti-fungal cream for athletes foot (as per NOK). Additional PMHX: unknown when diagnosed with Beta Thalassemia.	Levophed, Antihypertensives, NS, Propofol, NSBicarb, Fentanyl, Cardene, Zosyn, Levaquin, Labetalol, Vancomycin, Meperidine, Metoprolol, Clonidine, Norvasc, Hydralazine, Ancef, Keppra, Magnesium Sulfate, KCL, Heparin, Ativan, Morphine. Flushed and stored in UW solution.
HFC934	Female	54	C	CVA 2 nd ICH	Cigarettes- 2PPD X 15+ yrs. Marijuana (Smoked) Unknown amount; smoked daily X 40 yrs.	Emphysema, COPD X 1 yr. HTN- Diagnosed 6 mos. Prior; Compliant with meds. HTN medications, Metoclopramide, Proair HFA, Cymbalta, Lyrica, Levothyroxine, Amitriptyline, Premarin, Hydrocodone, Docusate Sodium, Calcium, Prenatal IPlus Tab, Simvastatin (as per NOK). PMHX: COPD, emphysema, back pain, fibromyalgia, bunion removal. Smoked marijuana and cigarettes daily. Pt was born with two uteruses, four ovaries, four kidneys (two of which were very small) and two livers. The extra organs were removed in 1979.	Fentanyl, Cisatracurium, Ancef, Albuterol, Hydrocortisone, Zyvox, Cefepime, Levothyroxine, Lasix, KCL, Furosemide, Zithromax, Rocephin, Propofol, Heparin, Magnesium Sulfate. Flushed and stored in HTK.
*HMC936	Male	60	C	Head Trauma 2 nd Blunt Injury	ETOH (Red Wine) 1 glass 4 X / wk X 30 yrs.	Enlarge Prostate med. (Unknown name of med), Multivitamins (as per NOK).	Propofol, Dobutamine, Neosyneprhine, Levophed, Insulin, Zosyn, Ancef, Vecuronium, Mannitol, Chlorhexidine, Chlorhexidine Gluconate, Docusate, Senna Syrup, Sodium Chloride, Phenylephrine, Acetaminophen, Calcium Gluconate, Fentanyl, Hydralazine, Hypromellose, Labetalol, Magnesium Sulfate, Morphine, Ondansetron, Ophthalmic Lubricant, Oxycodone, Potassium, Potassium Chloride. Flushed and stored in Custodial.

For Research Use Only. Not for use in diagnostic or therapeutic procedures.

© Discovery Life Sciences

Discovery Life Sciences
 6 Henshaw Street
 Woburn, MA 01801
 Tel: (866) 838-2798
info@dls.com
<https://www.dls.com/>

UltraPool™ HLM 150

Donor Information

Specimen	Gender	Age (years)	Race	Cause of Death	Social History	Medical History	Medication given during Hospitalization
*HMC937	Male	40	C	Anoxia 2 nd Cardiovascular	Cigarette: 1-1 1/2 PPW; Unknown duration' Chewing tobacco: Unknown amount, duration, frequency; Pt chewed more than smoked; Quit both 6 mos ago. ETOH (Beer) 2-3 beer- 12 pk at times, since high school; Quit less than 6 mos ago. Marijuana (Smoked) Occasional use off and on since High School, Unknown amount or frequency.	Stomach pain 8 yrs ago; Colonoscopy-Neg. Oxycodone, robaxin, Gabapentin, Zolof, Ibuprofen, Nerve medicine for back (as per NOK).	Epinephrine, Narcan, KPhos, Labetalol, KCl, MgSO ₄ , Fluconazole, Ancef, Vasopressin, Levophed, Neosynephrine, Levothyroxine, Solumedrol, Mannitol, Heparin. Flushed and stored in SPS-1
HMC938	Male	54	C	Anoxia 2 nd Cardiovascular	Non-smoker; Non-drinker	HTN; Diagnosed 3 yrs ago; Non-compliant w/ meds.	Vasopressin and epinephrine. Heparin, Nitroglycerin, Phenytoin, Zosyn, Sodium Phosphate, Propofol, Potassium Chloride, Fentanyl, Calcium Chloride, Cardizem, Lasix. Flushed and stored in HTK.
HMC941	Male	65	C	CVA 2 nd ICH	Cigarettes: 1/2 PPD for < 10 yrs; Quit in his 20's. ETOH (Beer, Wine) drank socially <3 drinks / wk; Unknown duration. Cocaine (not crack); THC; Unknown amount or duration; NOK states that it was a couple of time in college.	Seizures; During his late 20's-30's; Not taken meds for 30 yrs.	Cardizem, Baclofen, Vasopressin, Calcium Chloride, Labetalol, Potassium Chloride, Potassium Phosphate, Albuterol, Solu-Medrol, Insulin, D50, Levothyroxine, Naxolone, Zosyn, Sodium Bicarbonate, Dopamine, Levophed, Heparin, Mannitol. Flushed and stored in HTK.
HMC942	Male	56	C	CVA 2 nd ICH	Non-smoker; Non-drinker	HTN; Diagnosed 10-12 yrs ago. DM II Diabetes; Diagnosed 10 yrs ago; Had not taken meds in 2 yrs. Bowel obstruction in 2011. Losartan, Vitamin Z, Vitamin B, Vitamin D (as per NOK).	Levophed, Solumedrol, Insulin, Antihypertensives, DDAVP, Vasopressin, Cardene, Zosyn, Meropenem, Levaquin, Heparin. Flushed and stored in UW solution.
HMC943	Male	26	C	Anoxia 2 nd GSW	Chewing Tobacco: 1 bag / 2 wks X 4 yrs; Other Tobacco products: unknown name, frequency, and duration. ETOH (wine cooler) 1 wine cooler 1-2 x /mos X 2 yrs.	No meds taken at home (as per NOK).	Dopamine, Heparin. Flushed and stored in HTK.

For Research Use Only. Not for use in diagnostic or therapeutic procedures.

© Discovery Life Sciences

Discovery Life Sciences
 6 Henshaw Street
 Woburn, MA 01801
 Tel: (866) 838-2798
info@dls.com
<https://www.dls.com/>

UltraPool™ HLM 150

Donor Information

Specimen	Gender	Age (years)	Race	Cause of Death	Social History	Medical History	Medication given during Hospitalization
*HFC944	Female	44	C	CVA 2 nd ICH	Cigarettes- 1/2 PPD X 30 yrs. ETOH (Liquor) socially < 3 drinks; Unknown duration and frequency.	Asthma as a child. HTN X 5 yrs; Unknown if on medication or if compliant w/ meds. Tylenol PRN- nearly every day for migraines (as per NOK).	Albuterol, Insulin, Furosemide, Methylprednisone, Zosyn, Potassium Chloride, Duoneb, Solumedrol, Naloxone, Potassium Phosphate, Heparin, Mannitol. Flushed and stored in HTK.
HFC945	Female	52	C	Anoxia 2 nd Cardiovascular	Cigarettes- 1 PPD X 30 yrs. ETOH (Beer) Unknown amount, frequency, and duration although it is stated by NOK that it was socially or for special occasions only.	DMII Diabetes- Diagnosed at age 21; Controlled through PO meds and IV Insulin injections; Compliant w/ meds although glucose level was difficult to manage. (Names unknown to family) DM meds, Cholesterol meds, and Calcium supplement (as per NOK). PMHX- Seizures; Diagnosed w/ petite mal seizures between age 6 - puberty. During this time Pt had 2 seizures, was treated w/ medication, and outgrew them once she began puberty; Pt was told she had a narrowed blood vessel that apparently grew w/ age. No recurrence. Pt was not taking medication.	Propofol, Heparin, Morphine. Flushed and stored in HTK.
*HMC947	Male	68	C	CVA 2 nd ICH	Cigarettes: 1 PPD X 40 yrs; Quit in 2000. ETOH (Gin and Tonic) 1-2 drinks/ week X 48 yrs; Quit 1 yr ago.	Sleep Apnea w/ Cpap; Diagnosed in 2006; Unknown if compliant w/ Cpap. Massive MI in 2006 with Bypass. HTN; Diagnosed in 2007; Unknown if compliant w/ meds. DM Type I Diabetes; Diagnosed in 2007; Unknown if compliant w/ meds. Diverticulitis; Diagnosed in 2012 and treated w/ antibiotics in July and August; F/U colonoscopy 2 wks prior. Kidney Stones in 2012. ASA, Iron, Multivitamin, Januvia, Insulin, Clotrimazole-betamethasone, Omeprazole, Metoprolol Tartrate, Metformin, Simvastatin, Novolog, Lantus, Vitamin D, Depo-Testosterone (as per NOK).	Levophed, Insulin, T4, Heparin, Vecuronium, Decadron, Keppra, Levaquin, Protonix, Solumedrol, KPhos, CaCl. Flushed and stored in UW solution.
HMC949	Male	50	C	Anoxia 2 nd Cardiovascular	Cigarettes: 1/2 PPD since age 13; Quit 2 yrs ago.	Rheumatoid arthritis in knees and back. Macular degeneration. COPD x 4 yrs; Unknown if on med or compliant w/ meds. CAD x 5-6 yrs; Med controlled; Unknown if compliant w/ meds. HTN x 5-6 yrs; Med controlled; Unknown if compliant w/ meds. DM Type II Diabetes; Diagnosed in 2006; Medication controlled; Unknown if compliant w/ meds. Nicotine gum, Benadryl, Cyproheptadine, Vicodin, ASA, Valium, Divalproex, Gabapentin, PRimidone, Paxil, Quetiapine, metoprolol, Amlodipine, Isosorbide Mononitrate, Atorvastatin, Lasix, Insulin, Metformin, Albuterol, MDI (as per NOK). Pt took anti-malaria medication prophylactically for military.	Versed and Fentanyl drips. Levophed, Dopamine, T4, Insulin, Antihypertensives, Vasopressin, Heparin, Rocuronium. Additional PMHX- COPD, bipolar, PTSD, recurring leg infections, cyst removed from left wrist (non cancerous), wisdom teeth removed, HTN, DM, Angiogram following removal of 9 teeth, chronic ear infections. Flushed and stored in SPS-1.

For Research Use Only. Not for use in diagnostic or therapeutic procedures.

© Discovery Life Sciences

Discovery Life Sciences
6 Henshaw Street
Woburn, MA 01801
Tel: (866) 838-2798
info@dls.com
<https://www.dls.com/>

UltraPool™ HLM 150

Donor Information

Specimen	Gender	Age (years)	Race	Cause of Death	Social History	Medical History	Medication given during Hospitalization
*HMC950	Male	50	C	Head Trauma 2 nd Blunt Injury	Tobacco (Chew): x 10 yrs; Last time used 10 yrs ago; Unknown amount used or frequency. ETOH (beer) 6-12 cans/ week. Marijuana (Unknown route) Used recreationally; Unknown about used, frequency or duration.	Asthma; Diagnosed 3 yrs ago; unknown if compliant w/ meds. HTN x 15 yrs; Unknown if compliant w/ meds. GERD x 5 yrs; Bloody Stools occasionally, non diagnosis for it. Bystolic, Divalproex sodium, Atorvastatin, Ventolin HFA, Pulmicort Flexhaler, Omeprazole, Prednisone, Androgel, Fexofenadine HCL, Plavix, Tramadol, Nortriptyline (as per NOK).	KPhos, Levophed, Potassium, Magnesium, DDAVP, Zosyn, Fentanyl, Rocuronium, KPhos, Heparin. Flushed and stored in SPS-1.
HFC952	Female	40	C	Anoxia 2 nd Cardiovascular	Non-smoker, non-drinker. Marijuana X 5 mos from Sept. 2011- Jan 2012 for control of nausea; Unknown route, amount, and frequency.	Coronary Artery Bypass Graft in 2003; Quadruple Bypass; Unknown if compliant w/ meds. HTN X 10 yrs; Unknown if compliant w/ meds. Seizures; Diagnosed 2 yrs. Ago; Unknown if compliant w/ meds. DM Type I Diabetes; Diagnosed at age 7; Unknown if compliant w/ meds. Severe Gastroparesis for the past 10 yrs; Frequent vomiting and Diarrhea; Unknown duration and frequency. Kidney Failure; Diagnosed in 2008 d/t DM Type I; Pt was on Dialysis since 2008. Insulin, Lantus, Novolog, Levothyroxine, Bystolic, Simvastatin (as per NOK). PMHX; CAD, ESRD with dialysis three times per week, CVA, Hypothyroid, blindness in one eye d/t retinopathy to DM, secondary Hyperparathyroidism, Coronary Artherosclerosis, Hyperlipidemia, Esophagitis, Epilepsy, Jejunostomy feeding tube, Cesarean section in March 2012, Dialysis port right chest, eye surgery (right eye prosthetic).	Insulin, Levophed, Dopamine, Cardene, Labetalol, Solumedrol, Cefepime, Fluconazole, Levaquin, Protonix, Fentanyl, Lispro, Lantus, Lovenox, Zemplar, Heparin, Rocuronium. Flushed and stored in HTK.
HMC953	Male	76	C	CVA 2 nd ICH	Cigarettes: 1/2 PPD x 30 yrs; Quit in 1972; Occasional Cigars; Unknown amount, frequency, and duration. ETOH (Wine) <2 drinks / day x 40 yrs.	Quadruple Bypass in 1991. HTN x . 10 yrs; Compliant w/ meds. GERD; Unknown when diagnosed; Bloody stools resulting in a colonoscopy that revealed diverticulosis; Compliant w/ treatment. Kidney Stones; Incidence greater tan 10 yrs ago. PMHX; HTN, TIA, CABG, and had prostate removed in 2000 s/p cancer. Plavix, HCTZ, Ramipril, Lipitor, Xanax, Aspirin, Vitamins, Calcium, Chondroitin (as per NOK).	Neosynephrine, Solumedrol, T4, DDAVP, Ancef, Heparin, Zemuron. Flushed and stored in UW solution.
HFC955	Female	66	C	CVA 2 nd ICH	Cigarettes- 1 PPD every 2-3 days X 33 yrs. Quit 10 yrs ago. ETOH (Beer, Wine) rarely; Unknown amount, frequency, or duration	Multivitamin, Vitamin E, and Celexa (per NOK).	IV Fluids with Steroids, T4, Insulin, and Antihypertensives, Rocuronium, Metropolol, Cardene, and Vasopressin. Flushed and stored in UW solution.
HFC957	Female	36	C	CVA 2 ND to Stroke	cigarettes- 1/2-1 Pack per week for the last 4 yrs.	Hypertension- Diagnosed 1 yr. prior; pt compliant w/ meds. Lisinopril (as per NOK).	Heparin. Flushed and stored in UW solution.

For Research Use Only. Not for use in diagnostic or therapeutic procedures.

© Discovery Life Sciences

Discovery Life Sciences
6 Henshaw Street
Woburn, MA 01801
Tel: (866) 838-2798
info@dls.com
<https://www.dls.com/>

UltraPool™ HLM 150

Donor Information

Specimen	Gender	Age (years)	Race	Cause of Death	Social History	Medical History	Medication given during Hospitalization
*HMC960	Male	57	C	ICH 2 nd to CVA	Cigarettes: 1/2 PPD x 45 yrs. ETOH (Beer) rare.	PMHX; Glaucoma; Recent sinus infection x 1 wk; diverticulitis- repair of "hole" in colon 1996; arthritis in hands - per NOK unknown type	Vasopressin, Ancef. Liver flushed with SPS-1.
*HFC961	Female	37	C	Anoxia / Stroke	Non-smoker; Non-drinker	Heart Disease; Diagnosis name unknown to historian. Unknown names and types (as per NOK). History of pacemaker and several embolic strokes.	KCL, Lasix, Levothyroxine, Insulin, Narcan, Solumedrol, D50, Tobramycin, D5 1/2 NS, Artificial tears, Duoneb, Imipenem, NS, Diuretics, T4, Vasodilators and Antihypertensives. Flushed and stored in SPS-1.
*HMC962	Male	60	C	CVA 2 nd ICH	cigarettes: 1 PPD x 10 yrs; Quit 2-3 yrs ago. eTOH (Beer) 5-6 Beers daily x 10 yrs; Sober for 10 yrs but was a heavy drinker x 15 yrs. Marijuana (smoked) 1 time as a teenager; Unknown amount used.	HTN x 5-6 yrs; Non-compliant w/ meds. HCTZ, Amlodipine, Aspirin (as per NOK)	Dopamine, Neosynephrine, Soluortef, T4, Insulin, Heparin, Mannitol, Rocuronium, Dexamethasone. Flushed and stored in SPS-1.
HFC963	Female	74	C	CVA 2 ND to Stroke	1/4-1/2 PPD on and off; NOK unsure how long but Pt has not smoked in 5 yrs. 1/2 glass of wine during holiday visits to the UK annually X 54 yrs.	COPD diagnosed last year, Manic Depressive and Bipolar, compliant w/ Prozac. DM Type II diabetes; compliant with Insulin 4 x / day for 30 yrs. Pt. occasionally experienced persistent diarrhea 2 nd to diet. Kidney Stones when Gallbladder was removed 42 yrs. Ago. Prozac (as per Pt's chart). Hysterectomy 30 yrs prior, Cholecystectomy 42 yrs prior, and Pancreatectomy 30 yrs ago d/t yellow jaundice and Pancreatitis.	Vasopressin, Neosynephrine, Levophed, D50, Solumedrol, Lasix, Antihypertensives, Heparin, Mannitol, Vecuronium. Flushed and stored in SPS-1
*HFC964	Female	64	C	CVA 2 nd ICH	Cigarettes- 1 PPD X 20 yrs; Quit 24 yrs ago. ETOH (Wine) 3-4 nights/ wk on and off through adulthood; Unknown amounts but historian stated that she did not drink excessively.	Worked up for pulmonary nodule since 2007, PET scan and CT scans did not indicate malignancy; 7 mm LUL nodule representative of a benign calcified granuloma. Atrial Fib (required cardioversion), mitral valve murmur X 30 yrs; MVR w St. Jude's valve and AV pacer done in 2008. HTN; Unknown when diagnosed but Pt. was compliant w/ meds. Amoebic dysentery x 2 while she was in the Peace Corps in Guatemala; Unknown when diagnosed. Metoprolol, Quinapril, Coumadin, HCTZ, ASA, Simvastatin, MVI, Vitamin C (as per NOK). Hysterectomy > 25 yrs ago.	Insulin, Potassium Chloride, Clindamycin, Maxipime, Lasix, Decadron, Levophed, MagSulfate, Insulin, Vecuronium, Neosynephrine, Heparin. Flushed and stored in HTK.
HMC965	Male	46	C	Anoxia 2 nd to Asphyxiation	cigarettes: 3 PPD x 20 yrs. ETOH (Beers) 12 beers/ day x 18 yrs. Marijuana (Smoke) Occasionally 1-2 x/ yr x2 yrs; Last time was 1982.	Anxiety w/ stress causing chest pain; Unknown diagnosis or when it began. Multivitamin, vitamin D (as per NOK).	Epinephrine, Methylprednisolone, Hydrocortisone, Insulin, Vasopressin, Sodium Phosphate, Heparin. Flushed and stored in UW solution.

For Research Use Only. Not for use in diagnostic or therapeutic procedures.

© Discovery Life Sciences

Discovery Life Sciences
6 Henshaw Street
Woburn, MA 01801
Tel: (866) 838-2798
info@dls.com
<https://www.dls.com/>

UltraPool™ HLM 150

Donor Information

Specimen	Gender	Age (years)	Race	Cause of Death	Social History	Medical History	Medication given during Hospitalization
*HFC966	Female	65	C	CVA 2 nd ICH	3 PPD X 49 yrs; Quit on 11/19/2012. ETOH (Beer) 6 pk / day x >40 yrs.	Asthma; Unknown when diagnosed or if compliant w/ meds. HTN x > 10 yrs; Unknown if compliant w/ meds. UTI 1 yr ago. Terazosin, Omeprazole, Loratidine, Fluticasone, Pro-Air Inhaler, CoQ-10, Vitamin C, Omega 3, Absorbable Calcium, Biotin, Alfalfa, B12, Potassium Citrate (as per NOK).	Methylprednisone, Hydrocortisone, Insulin, Vasopressin, Dopamine, Invanz, KCl. Flushed and stored in UW solution.
*HFC969	Female	59	C	HT 2 nd SIGSW	Cigarettes- 1 1/2 PPD X 40 yrs. ETOH (unknown type) 2 drinks / week; unknown duration.	Asthma; Unknown when diagnosed or if compliant w/ meds; COPD; Unknown when diagnosed or if compliant w/ treatment. HTN; Unknown when diagnosed; Compliant w/ meds. Neuropathy; Unknown when diagnosed or if compliant w/ meds. Additional PMHX: Venous Insufficiency and Thromboangitis Obliterans.	Vancomycin, Lasix, Levophed, Lasix, Ancef, Vancomycin, Heparin. Flushed and stored in SPS-1.
*HFC970	Female	75	C	CVA 2 nd ICH	Cigarettes- Unknown amount or frequency X 30 yrs; Quit 20 yrs. ago. ETOH (Wine, Liquor) 1 glass of wine or hard liquor now and again; Unknown frequency or duration.	Asthma; Diagnosed as a child; COPD; Diagnosed 3 yrs. prior; Unknown if compliant w/ meds. Kidney Stones; Diagnosed 5 yrs. Ago and were treated w/ laser therapy. Spiriva (as per NOK).	Neosynephrine, Dopamine, T4, Vasopressin, Cipro, Heparin. Flushed and stored in HTK.
HMC972	Male	56	C	Anoxia	cigarettes: 2 PPD x 45 yrs. ETOH (Beer) Occasionally drank since Pt was teenager; Unknown amount or frequency.	Emphysema; Unknown when diagnosed or if compliant w/ treatment. Schizophrenia; Unknown when diagnosed or if compliant w/ treatment. Med for Schizophrenia (as per NOK).	Acetaminophen, Albuterol HFA, Amiodarone, Atropine, CaGluconate, Ceftriaxone, Chlorhexidine, DDAVP, Diphenhydramine, Dopamine, Doxycycline, Lasix, Hydralazine, Insulin, Labetalol, Solumedrol, Versed, Morphine, Levophed, Phosphate Sodium, Phosphate Potassium, Potassium Chloride, Vasopressin. Flushed and stored in SPS-1.
HMC974	Male	33	C	HT 2 nd Blunt Injury	Cigarettes: 1 PPD x 4 yrs. ETOH (Beer, Liquor) drank daily; Unknown amount or duration.	No meds taken at home (as per NOK).	Sodium Bicarbonate, Furosemide, Mannitol, Albuterol, Potassium Phosphate, Furosemide, Zosyn, Levophed, Propofol, Fentanyl, Heparin. Flushed and stored in HTK.
*HMC975	Male	26	C	HT 2 nd Blunt Injury	Dipped Skoal; 2-3 x / wk x 7 yrs. ETOH (Beer) 2-3 at a time, 2-3 x / week x 5 yrs.	No meds taken at home (as per NOK).	KCl, Neosynephrine, T4, Ancef, Solucortef, Zosyn, Lasix, Vancomycin, Albuterol, Keppra, Levophed, Mannitol, Morphine, Fentanyl, Ativan, Rocuronium, Albumin. Flushed and stored in SPS-1.
*HFC979	Female	27	C	Head Trauma 2 nd Blunt Injury	Cigarettes- 1 PPD X 5 yrs. Wine- Occasional drinker < 2 drinks per day (as per NOK).	Donor did not take any medications (as per NOK).	Vasopressin, Levophed, Ancef, FFP, D5W. Flushed and stored in HTK.

For Research Use Only. Not for use in diagnostic or therapeutic procedures.

© Discovery Life Sciences

Discovery Life Sciences
6 Henshaw Street
Woburn, MA 01801
Tel: (866) 838-2798
info@dls.com
<https://www.dls.com/>

UltraPool™ HLM 150

Donor Information

Specimen	Gender	Age (years)	Race	Cause of Death	Social History	Medical History	Medication given during Hospitalization
*HFC982	Female	62	C	CVA 2 nd ICH	Cigarettes- 1 PPD X 20 yrs.	Pneumonia; Recently Diagnosed w/ community acquired Pneumonia; Unknown if compliant w/ meds. CHF and Afib x 30 yrs; Pacemaker inserted in 2000; Compliant w/ meds. HTN x 30 yrs; Compliant w/ meds. Memory loss with in the last 2 mos with headaches. Cervical Cancer; Diagnosed 25 yrs ago; Last check up was 5 mos ago and no treatment was required. Coumadin, Heart Meds, Furosemide, Morphine as per NOK). Chronic back pain; Thoracic and Lumbar fusions.	Levofloxacin, Levophed, Pitressin, KPhos, Milrinone, Heparin. Flushed and stored in UW solution.
*HMC984	Male	70	C	HT 2 ND Blunt Injury	Cigarettes: 2 PPD x 20 yrs; quit 40 yrs ago (NOK). ETOH (Beer) 6 pk daily; stopped 6 yrs ago (NOK). Marijuana; unknown amounts from 1967-1969 (NOK).	HTN x 5 yrs; med compliant. NIDDM x 1 yr. Kidney Stones. Losartan, Terazosin, Metformin, Prilosec, Gabapentin, Aspirin, Iron, Loratadine, and Vitamins on a regular basis (per NOK).	PRBCs, D5, Vasopressin, Levophed, KCL, Bumex, Solumedrol, Levothyroxine, Ancef, T4, Dopamine, Insulin, Albuterol, Keppra, Nembutal, Magnesium Sulfate, Cefepime, Vancomycin, Docusate, Famotidine, Versed, Etomidate, Vecuronium and Platelets. Flushed and stored in SPS-1
*HFC987	Female	63	C	CVA 2 nd ICH	Cigarettes- 5-10 / day x 40 yrs; Pt did cut down within the last 2 yrs although amount unknown. ETOH (Brandy) 3-4 drinks / day x 10 yrs.	Afib; Unknown when diagnosed or if compliant w/ treatment. HTN x 5 yrs; Unknown if Pt was compliant w/ meds. Nifedipine, MVI, Amiodarone, Zetia, MgOxide, Lotensin, Synthroid, Coreg, Coumadin, Aspirin (as per Pt's chart)	Labetalol, Colace, MgHydroxide, Dulcolax, Novolog, D50W, Artificial Tears, Protonix, Combivent, Cardene, KCl, Zofran. Flushed and stored in UW solution.
HMC988	Male	52	C	Anoxia 2 nd Cardiovascular	Cigarettes: 1 PPD x 40 yrs; Quit 2-3 weeks ago. ETOH: "black velvet" 1 pint daily x 15 yrs; Quit 13 yrs ago.	mild COPD; Diagnosed 2 yrs ago. Sleep apnea; wears a bi-pap at night x 15 yrs. HTN x 5 yrs; med compliant. Diabetes x 2-3 yrs; med compliant. Baclofen, Norco, Claritin, Glucophage, Methadone, Inderal and Imitrex on a regular/daily basis (as per NOK).	Levophed, Vasopressin, Solumedrol, Lasix, T4, Insulin, Antihypertensives, Heparin, KPhos, Magnesium Sulfate, KCL, Acetylcysteine, Calcium Gluconate, Sodium Bicarbonate, Labetolol, Clindamycin, D50, Narcan, Insulin, Vecuronium, Unasyn, Potassium Chloride, Hydralazine, NS, Atrovent and Mucomyst. Flushed and stored in HTK.
HMC992	Male	30	C	HT 2 nd Blunt Injury	ETOH (Liquor, Beer) Fruity Drinks; Maybe 1 x / mos but very rare; Beer; Unknown amount, frequency, or duration.	DM Type II; Diagnosed in 2008 x 1 yr; Pt lost weight and it resolved itself. No meds taken at home per NOK.	Albuterol, KCl, K-Phos, Zosyn, Vancomycin, Insulin, Fentanyl, Lovenox, Ancef, Nicardipine, Levophed. Flushed and stored in HTK.

For Research Use Only. Not for use in diagnostic or therapeutic procedures.

© Discovery Life Sciences

Discovery Life Sciences
6 Henshaw Street
Woburn, MA 01801
Tel: (866) 838-2798
info@dls.com
<https://www.dls.com/>

UltraPool™ HLM 150

Donor Information

Specimen	Gender	Age (years)	Race	Cause of Death	Social History	Medical History	Medication given during Hospitalization
*HFC993	Female	65	C	Anoxia 2 nd Cardiovascular	Cigarettes- 2 PPD x 45 yrs. ETOH (Beer, Liquor) a 5 th of Scotch daily from teen yrs to age 40; Hasn't had alcohol x 20 yrs. Heroin (IV) over 15 yrs ago; Unknown frequency.	COPD; Unknown when diagnosed or if compliant w/ treatment; Emphysema; Unknown when diagnosed or if compliant w/ treatment; Pneumonia. 2 CVA; Unknown when diagnosed or if compliant w/ treatment; Unknown date of when Pt underwent Carotid Endarterectomy; MI; Unknown when diagnosed or if compliant w/ treatment; Pt was admitted for chest pain and had her Right coronary Artery stented 5 days prior to last admission. HTN x 20 yrs; Unknown if compliant w/ treatment. Seizures; Unknown when diagnosed or if compliant w/ treatment (as per Pt's chart). DM Type II Diabetes x 25 yrs; Unknown if compliant w/ treatment. GERD; Unknown when diagnosed or if compliant w/ treatment. Advair, Albuterol, Aldactone, Bactrim DS, Lisinopril, Lortab, Lyrica, Metformin, Mucinex, Nitrostat, Paxil, Plavix, Prilosec, Senokot, Spiriva, Medication for Osteoporosis- name unknown to Historian (as per NOK).	Dopamine, Dobutamine, Solumedrol, Mannitol, T4, Insulin, Vasodilators, Vasopressin, Heparin, Rocuronium, Dexamethasone. Flushed and stored in HTK.
HMC996	Male	70	C	HT 2 nd SIGSW	Cigarettes: 1/2 PPD x 4-5 yrs; Cigar: 2-3 / day x 20 yrs. Marijuana (Unknown route) x 1, 2-3 yrs ago.	Pneumonia; Diagnosed 15 yrs ago. Multiple Sclerosis x 30 yrs; Progressive loss of gait d/t MS. Constipation & Diarrhea x 2-4 yrs w/ incontinence. Oxycodone, Fentanyl Patches, Lidocaine Patch, Sleeping Pills, Vitamin C, D, E, Fish Oil, MVI (as per NOK).	Epinephrine, Levophed, T4, Solumedrol, Insulin, Neosynephrine, Hydralazine, Lasix, KCl, CaGluconate, Soluortef, Cefazolin, MgSulfate, Propranolol, DDAVP, Dilantin, D5W, Pepcid, Vecuronium, Midazolam, Fentanyl, Succinylcholine, Etomidate, Senna, Heparin. Flushed and stored in SPS-1.
*HFC997	Female	48	C	CVA 2 nd ICH	Cigarettes- 1/2 PPD X 28 yrs. ETOH (Mixed Drinks) Unknown amount, frequency; Pt consumed in her 30's, would increase with Mania and subside when depressed/normal. Cocaine in her 20's; Unknown route, amount, frequency; Speed (Snorted) in her 20's; Unknown amount and frequency; Pills in her 20's; Unknown amount or frequency.	HTN; Diagnosed in 2011; Compliant w/ treatment. Dementia; Diagnosed in 2011; Unknown if compliant w/ meds; Depression; Unknown when diagnosed or if compliant w/ treatment; Bipolar Disorder; Unknown when diagnosed or if compliant w/ treatment. MVI, Lithium, Seroquel, Calcium Supplements, Thyroid Medication, HTN Medication, Klonopin, Docusate, T4, Valacyclovir, Venlafaxine, Risperidone, Simvastatin (as per Pt's chart). Additional PMHx: HLD, hypothyroidism.	Dopamine, Levophed, Famotidine, Keppra, Vasopressin, Dextrose, T4, Insulin, DDAVP, Heparin. Flushed and stored in HTK.
HFC999	Female	73	C	CVA 2 nd ICH	Non-smoker; Non-drinker	OTF "Bone Pill" from Health Food Store (as per NOK).	Levophed, Vecuronium, Etomidate, Versed, Fentanyl, Vasopressin, Dopamine, Solumedrol, T4, Insulin, DDAVP, Cipro, Heparin. Flushed and stored in HTK.

For Research Use Only. Not for use in diagnostic or therapeutic procedures.

© Discovery Life Sciences

Discovery Life Sciences
6 Henshaw Street
Woburn, MA 01801
Tel: (866) 838-2798
info@dls.com
<https://www.dls.com/>

UltraPool™ HLM 150

Hazard Warning:

- This microsome preparation was prepared from freshly frozen human tissues. All donor tissues have tested negative for pathogens by PCR for the following: HIV I/II, HTLV I/II, CMV, HBV and HCV, however, we recommend that this material be considered a potential biohazard.
- Donors with positive serology for CMV are identified in the donor demographic sheet with a single asterisk. Donors with CMV serology unknown are identified with a double asterisk. Donors CMV negative for serology are unmarked.

Safety Information:

- Chemical hazard assessment indicates this product is not hazardous, therefore no SDS (Safety Data Sheet) is provided.
- This product was prepared from freshly frozen human tissues. All donor tissues have tested negative for pathogens by PCR for the following: HIV I/II, HTLV I/II, HBV and HCV. Donors with positive serology for CMV are identified in the donor demographic sheet with a single asterisk. Donors with CMV serology unknown are identified with a double asterisk. Donors CMV negative for serology are unmarked.
- All persons handling human-derived materials must use Universal Precautions in accordance with the US OSHA Bloodborne Pathogens Final Standard and/or the European Council Directive on the protection of workers from risks related to exposure to biological agents at work (90/679/EEC) and it's relevant European National Transpositions.



20 June 2024

Quality Assurance

Date

For Research Use Only. Not for use in diagnostic or therapeutic procedures.

© Discovery Life Sciences