

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.	38301
Qty./Package	0.5 mL
Protein Content	20 mg/mL in 250 mM sucrose
Storage Conditions	Store at -80°C
Release Date	2024 September
Expiration Date	2034 May

Assay Results

Enzyme Measured	Assay	Enzyme Activity [in pmol/(mg x min)]
Total P450	Omura and Sato	330 pmol/mg
OR	Cytochrome c Reductase	350 nmol/(mg x min)
Cytochrome b ₅	Spectrophotometric	400 pmol/mg
CYP1A2	Phenacetin O-deethylase	590
CYP2A6	Coumarin 7-hydroxylase	1000
CYP2B6	Bupropion hydroxylase	270
CYP2C8	Paclitaxel 6 α -hydroxylase	280
CYP2C9	Diclofenac 4'-hydroxylase	2600
CYP2C19	(S)-Mephenytoin 4'-hydroxylase	60
CYP2D6	Bufuralol 1'-hydroxylase (The amount of activity inhibited by 1 μ M quinidine.)	90
CYP2E1	Chlorzoxazone 6-hydroxylase	2300
CYP3A4	Testosterone 6 β -hydroxylase	5700
CYP4A11	Lauric acid 12-hydroxylase	1200
FMO	Methyl p-Tolyl Sulfide Oxidase	1900
UGT1A1	Estradiol 3-Glucuronidation	1000
UGT1A4	Trifluoperazine N- Glucuronidation	680
UGT1A6	Serotonin Glucuronidation	15000
UGT1A9	Propofol Glucuronidation	3400
UGT2B7	AZT Glucuronidation	1800
Western Blot Results		
CYP3A4	Western Blot	110 pmol/mg
CYP3A5	Western Blot	26 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.

© DLS

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
HFC1	Female	31	C	MVA	Tobacco use; 1 PPD x 15 yrs.	No medications	Not available.
*HMC32	Male	20	C	MVA	Marijuana use.	Unknown	Dilantin, Morphine, Ancef.
HFC56	Female	57	C	Anoxia	Not available.	Amoxicillin, HTN, Pulmonary Fibrosis, Arthritis, Diabetes, Insulin Injections	Dopamine
HFC82	Female	61	C	Intracranial Bleed	Not available.	Hypertension; non-compliant w/ meds.	Not available
HFC101	Female	54	C	MVA	Not available.	Healthy Normal	Not available.
HMC121	Male	59	C	GSW	Social Alcohol. Tobacco use; 2 PPD x 40 yrs.	Skin cancer, basal cell, unconfirmed lung cancer, Elavil, Lithium.	Dopamine, Levophed, Atropine, Epinephrine, Lasix, DDAVP, Lidocaine, Ancef.
HMC181	Male	50	C	Head trauma, MVA	Alcohol use: holidays. Marijuana: age 18.	Maalox, kidney stones-MV, Prolapse.	Neosynephrine, Vasopressin, Fentanyl, Insulin, Pentobarb, Gentamycin, Ancef, Vancomycin.
HMC184	Male	21	C	Respiratory failure	Smoked: 1 PPD x 2 yrs. 5-6 12-oz. beers/week, jail for one night for DUI	Adrenoleukodystrophy, radiation, chemotherapy + stem cell transplant /2006. Kidney failure and dialysis during surgery, took dexamethazone, Florinef, Paxil, and Resperidol, diffuse alveolar damage.	Crystalloids, colloids, blood products, hydromophone, furosemide, calcium.
HFC189	Female	54	C	CVA	Alcohol once or twice a year.	MVI and Calcium.	Nipride, Insulin, Norepinephrine, Vasopressin, Unasyn, Levaquin, Ancef, Heparin, Levophed
*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, Vancomycin, Etomidate, Lidocaine, Tetanus, Vecuronium.
*HFC203	Female	46	C	CVA	Smoked Metamphetamine x 10 yrs	Benign	NS, T4, Insulin, Dopamine, Neosynephrine, K+Acetate, Albuterol, DDAVP, Ancef, D5W, KPhos and LR

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

© DLS

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
*HFC217	Female	45	C	Anoxia	Non-smoker, Non-drinker.	Cognitive Disorder; moderate mental retardation since birth. Omeprazole, Mucinex, Ibuprofen, Colace, Acetaminophen, Albuterol, and MOM (as per NOK).	Heparin, Regitine, Omeprazole, Mucinex, Bisacodyl, Ibuprofen, Docusate Sodium, Adcetaminoph3en and Albuterol. Flushed and stored in UW solution.
*HFC232	Female	66	C	CVA / Stroke	Non-smoker, Non-drinker.	Hypertension; Diagnosed 10 yrs ago. Nexium and anti-hypertension meds (as per NOK).	Dopamine, Tylenol, D5LR, and Narcan.
HFC237	Female	42	C	CVA	Non-smoker; Non-drinker	Cardiac Catheter one year ago, Distal LAD 80% greater than 50% RCA. Pulmonary Hypertension; compliant with Lisinopril and Catapress. Type I Diabetes; diagnosed at age 13; compliant with Novalin, NPH, and Regular Insulin,. ESRD; dialysis start one yr ago 3 q wk. Coumadin, LIsinopril, Catapres, Metoprolol, Reglan, Pepcid, Insulin, Lorazepam, Darvocet, and Tylenol (as per NOK).	Dopamine, Levophed, Neosynephrine, Insulin, Albuterol, Atrovent, D50, Heparin, KCL, Magnesium Sulfate, Narcan, Norcuron, Potassium Chloride, Protonix, Solumedrol, T4, and Vancomycin.
*HFC239	Female	73	C	CVA	Non-smoker; Non-drinker	Triple Bypass Surgery one yr prior. Diabetes; Diagnosed at age 50; treated with insulin.	Dopamine, Insulin, Vasopressin, Zosyn, KCL, Hydralazine, Solumedrol, Magnesium, Dopamine, Insulin Lasix, D50, NaHcO3, Potassium, DDAVP. Kphos, and Chloride. Flushed and stored in HTK.
HFC252	Female	66	C	CVA	Beer- very infrequently.	Pt had (+) TB test in 1960's, not tested since. CAD-Heart Stent placed. Hypertension; Diagnosed 40 yrs ago. Diabetes; Diagnosed 20 yrs ago; compliant with Glyberide. Gastroparesis and Diverticulitis. Multivitamin, Toprol, Glyburide, Lortab, Lisinopril, Reglan, Oxybutinine, ASA, Nexium, and Phenergan (as per NOK).	Heparin, Levophed, Dopamine, Neosynephrine and D51/2 NS. Flushed and stored in HTK.
HFC253	Female	63	C	CVA	Cigarettes; 30 yrs; Quit 5 yrs ago. Wine- moderate q d.	Hypertension. Plavix, Darvon, Zanaflex, Zolof, Lisinopril and Tylenol (as per NOK).	Neosynephrine, Pitressin, and Solumedrol.
HMC258	Male	42	C	Anoxia	Cigarettes; 1 PPD x 24 yrs. Liquor & Beer 2-3 drinks per month. Cocaine; 1 or 2 times when 21. Marijuana; quit 20 yrs ago.	Epiglottitis diagnosed 5 yrs ago. HBP diagnosed 5 yrs ago. Controlled with Diovan. Diovan, Viagra, Prednisone, Amoxycillin, Inhaler, Simvastatin (as per NOK).	Heparin. Flushed and stored in HTK.
*HFC259	Female	72	C	Head Trauma- Blunt Injury	Cigarettes; 1 PPD x 30 yrs.	Hypertension; Diagnosed 30 yrs ago. Controlled with oral medication. BP med unknown (as per NOK).	Insulin, Hydrocortisone, KCL, Vitamin K, Kkondremul, Sennakot, Pepcid, Protamine, Labetalol, Hydralazine, Heparin, Neosynephrine, Vecuromium. Flushed and stored in HTK.

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

© DLS

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
HMC263	Male	35	C	ICH	Beer; 2-3 twice monthly. Cocaine; Snorted less than once q month; last used 4 weeks ago. Marijuana; quit 5 yrs ago.	Asthma; Since childhood; No intubations. Flovent and Albuterol (as per NOK).	Dopamine, Solumedrol, Recombinant, Ancef, Mg Sulfate, Kphos, CaCl, Lasix, Solumedrol, Zosyn, D5NS, and Vasopressin.
*HMC269	Male	62	C	Head Trauma-MVA	Chewing Tobacco-Quit 5 yrs ago. Beer and Shots. Heavily-Quit 16 yrs ago.	Black Lung, HBP-Diagnosed 3-4 yrs ago. No neurological disease but pt has experienced some recent memory loss- starting 1 wk prior. Borderline Diabetes-Diet Controlled. GERD. Spironolactone, Cyanocobalamin, Benzonatate, Digoxin, Iron, Bupropion, Theravite, Bumex, Lopressor, Viagra, Memantine, Citalopram, Omeprazole, Donepezil, and levetiracetam (as per NOK).	Dopamine, Levophed, Insulin, Keppra, Crystalloids, Heparin, Ancef and Methylprednisolone. Flushed with UW.
*HMC270	Male	48	C	SIGSW	Cigarettes; 1 PPD x 30 yrs.	Meniere Disease for the last 12 years, syncope. Depression. Diarrhea, Stomach Cramps related to nerves. Vitamins, Fish oil, Syboxin, Cymalta, and Xanax (as per NOK).	Vasopressin, Insulin, Pepcid, Dextrose, Zofran, Ondansetron, Dopamine, Fentanyl, Norepinephrine, Sodium Chloride, Propofol, Calcium Chloride, Levothyrosine, T4, Pitressin, PRBCs and Magnesium Sulfate. Flushed and stored in UW solution.
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, Neosynephrine, Dopamine, Vasopressin, Kphos, Ancef, solumedrol, and Cleocin. Flushed with UW solution.
HMC282	Male	41	C	ICH-Stroke, MVA	Tobacco- 1/2 PPD X 25 yrs. Beer- 1/ mo since age 21. Cocaine/Heroin-Snorted Unsure of amount from 1980's to present.	Normal healthy	Dobutamine, Cacl, Insulin, Kphos, Levothyroxine, D50, Vasopressin, Solumedrol, Ancef, Mannitol, Levophed, Dopamine, Tylenol, Cerebyx, Dilaudid, PRBCs, Albumin 5%, 1/4 NS, 40 KCL, Heparin and Tridil. Flushed and stored in Viaspan.

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

© DLS

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
*HMC287	Male	76	C	ICH-Stroke	cigarettes- 1 1/2 PPD X 40 yrs. Quit 6 1/2 yrs ago. Scotch 2-3/week X 40 yrs. Quit February 2008.	Arrhythmia. Pace maker place 2/08. Congestive Heart Failure 12/06. Hypertension; Diagnosed 2 yrs prior. Dementia; Undiagnosed. Family attributes loss of memory due to head trauma throughout life as football player. Polyps; Diagnosed mid 80's. At least 50 polyps removed. Surgery in 1995 for removal of large polyp. Recurring Gout 1 1/2 yrs ago. BP meds, Adema meds, Zolof, and CHF meds (as per NOK).	Acetaminophen, Acetylcysteine, Calcium Gluconate, Dextrose, digoxin, Epinephrine, Esmolol, Furosemide, Heparin, Hydralazine, Levetiracetam, Mannitol, Metoprolol, Neosynephrine, Piperacillin, Vancomycin, NS, Levophed, Solu-Medrol, Zosyn, Albuterol, Insulin, Lasix and Vecuonium. Flushed and stored in Custodial.
*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
HFC308	Female	61	C	ICH-Stroke	Non-smoker; Non-drinker	Mild Seasonal Asthma. Diagnosed with Hypertension 30 yrs prior to death. Darvon, Ativan, Verapamil SR, Naproxen, and Detrol LA (as per NOK)	Neosynephrine, Vasopressin, Ancef, Fortaz, Heparin, Mannitol, and Lasix. Flushed and stored in UW solution.
*HFC316	Female	70	C	ICH-Stroke	Cigarettes: 1/2- 1 PPD X 53 yrs.	Carotid Artery surgery- 6 months prior to death. Diagnosed with hypertension 6 months prior to death. Alzheimer's: Diagnosed 4-5 yrs. Prior to death. Detached Kidney-Diagnosed 10 yrs. Prior to death. Namenda, Plavix, Crestor, Exelon, and Oxybutynin (as per NOK)	Levophed, 1/2 NS, Neosynephrine, Levothyroxine, K-Phos, Insulin, D50, Solumedrol, Ancef, Dopamine, .45 NS, 0.9% NACL, Albumin 5%, Crystalloids, Heparin, Vasopressin and Epinephrine. Flushed and stored in UW solution
*HFC329	Female	30	C	Anoxia-2 nd to Respiratory Failure	Non-smoker; Non-drinker	Rare interstitial lung disease; treatment started 10 yrs prior. Pt received bilateral lung transplant 2 yrs prior. Hypertension; diagnosed within the last 3-4 months; took HCTZ on and off. Hx of pulmonary HTN. Tacrolimus, Prednisone, Posaconazole, Bactrim, Topamax, Acyclovir, Allegra, Protonix, Trazodone, HCTZ, Azithromycin, Levaquin, Nasonex, Actenol, Advair, Xopenex, Tylenol, Frovatriptan, Calcium +D, Multivitamins and K-Dur (as per NOK).	Crystalloids, Insulin, Acryclovir, Azithromycin, Fexofenadine, Heparin, Hydrochlorothiazide, Dopamine, Lasix, Magnesium Sulfate, metolazone, Metronidazole, Potassium Chloride, Prednisone, Sertraline, Trazadone, Vancomycin and Zosyn. Flushed and stored in HTK.
*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.

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

© DLS

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
*HFC346	Female	52	C	ICH-Stroke	Cigarettes: 2 PPD X 15 yrs.	Hypertension diagnosed in 1992. Diabetes diagnosed in 1999. Insulin dependent X 4 yrs. Spironolactone, Nexium, Combivent, Crestor, Cardizem, CD-240 and Hydrocodone (as per NOK)	Insulin, Dopamine, Neosynephrine, Levaquin, Lasix, Solumedrol, Ancef, KCL, Heparin, Mannitol and Pavulon. Flushed and stored in UW solution.
*HFC351	Female	66	C	ICH-Stroke	Non-smoker; Non-drinker	Asthma/COPD- Unknown time/period. Cardiac disease- unknown time/period. Hypertension- unknown time/period. Diabetes- diagnosed 10 yrs prior. Insulin dependent X 6 yrs. Coumadin, Levothyroxine, Crestor, Plavix, Nexium, Cymbalta, Humulin, and Insulin (as per NOK)	1/2 NS, Labetalol, Levothyroxine, Insulin, Solumedrol, Esmolol, Nipride, 1/4 NS, Vecuronium, Levaquin, D50, Vitamin K, Heparin, and Crystalloids. Flushed and stored in UW solution.
*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, Phenergam, 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.
HMC383	Male	53	C	Head Trauma- SIGSW	Cigarettes: 1/2 PPD X 23 yrs. Quit in 2007. Beer/Wine: Socially.	Diabetes; Diagnosed 2 yrs ago. Tx with Glucophage. BP Meds, Vitamin C & E, Coumadin and Hydrochlorizide (as per NOK).	Crystalloids, Colloids, Neosynephrine, Vasopressin, Dobutamine, Levothyroxine, Insulin, Levophed.
HMC384	Male	64	C	ICH-Stroke	Cigarettes: 2 1/2 PPD X 50 yrs.	Hypertension; Diagnosed 15 yrs ago. Type II Diabetes; Diagnosed 2001. Problems urination at times. Advil and Aspirin (as per NOK).	Hespan, 1/2 NS, Heparin, Mannitol, Lasix, Epinephrine, Levothyroxine, Solumedrol, Electrolytes, Vasopressin and Cefazolin. Flushed and stored in UW solution.
*HFC408	Female	42	C	CVA	Cigarettes- 1-1.5 PPD X 28 yrs.	Hypertension- Diagnosed 1.5 yrs prior. Controlled w/ diet. Type II Diabetes- Diagnosed 1.5 yrs prior. Compliant with Metformin. Metformin, Ibuprofen, and Cholesterol meds (as per NOK)	Levothyroxine, Dopamine, Solucortef, Insulin, KCL, NaHCO, Calcium Gluconate, DDAVP, Epinephrine, NS, Vasopressin, KCL, MagSulfate, Solumedrol and Ancef. Flushed and stored in HTK.

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

© DLS

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
*HFC432	Female	40	C	ICH-Stroke	Non-smoker; Non-drinker	Pulmonary Embolism- Unknown when diagnosed. COPD- Unknown when diagnosed. Asthma- Diagnosed at age 17. Diabetes controlled with oral meds from 1996 - 2004. Resolved post Gastric Bypass. GI bleed X 1 yr status post gastric bypass 6/04. Coumadin, Oxycontin, Alprazolam, Promethazine, Lasix, Levothyroxine, Citalopram Hydrobromide, Potassium Chloride, Prenatal Vitamins, Fluticasone-Salmeterol inhaler and Albuterol inhaler (as per NOK)	1/4 NS with 20 KCL, D5 1/2 NS, D5W, LR, FFP/Plasma, Platelets, Dopamine, Neosynephrine, Ancef, Cardizem, Cerebyx, Desmopressin, Dextrose, Diltiazem, Esmolol, Fentanyl, Fosphenytoin, Hydralazine, Insulin, KCL, Labetalol, Lasix, Levothyroxine, Lorazepam, Magnesium Sulfate, Mannitol, Naloxone, Phenylephrine, Phenytoin, Solu-Medrol, T4, Vitamin K, Zosyn and Heparin, Flushed with LR, Additives- Solumedrol, Lidocaine, and Verapamil.
*HFC433	Female	82	C	Blunt Head Trauma	Non-smoker; Non-drinker	Atrial Fibrillation- Unknown when diagnosed. Hypertension- Diagnosed 7 yrs ago.	Neosynephrine, T4, Levothyroxine, Ancef, Crystalloids, Colloids, 5 % Albumin, Platelets, Heparin, Mannitol, and Lasix. Flushed and stored in UW solution.
HFC446	Female	76	C	ICH-Stroke	Cigarettes- 1-1.5 PPD X 62 yrs. Wine- socially.	Emphysema- Possible, but was never followed up on. ASA and Tylenol (as per NOK).	Levophed, Zosyn, D5 1/2 NS, D5W, NS, KCL, KPhos, Calcium Gluconate, Nicardipine, Heparin, Vancuronium, Insulin, LR and Vasopressin. Flushed and stored in HTK.
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.
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, Cefapime, Vancomycin, Potassium, Phosphate, Glycolpyrrolate, Insulin, Hetastarch, Heparin, and Levophed. Flushed an stored in HTK.

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

© DLS

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
*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.
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.
*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.

© DLS

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.
HMC518	Male	47	C	Single gun shot wound to head	1 PPD smoker, light drinker couple beers/wk. Overnight jail stays in past. Marijuana user in the past.	Healthy male	Dopamine, NEO, Pavulon, Heparin
HMC519	Male	70	C	Brain stem infarct	Normal.	Type II DM, controlled with oral insulin. Parkinson's. Appendectomy. Renal angioplasty 1990. Below knee (right) amputation due to emboli 1990. Breast lump. Degenerative joint disease, managed with non-steroidal anti-inflammatories. Pyloric ulcer, recent, duodenal ulcer. Recently hospitalized for fever, discharged 6/6/94. Vasotec, Prilosec, Symmetrel, Dulcolax, Carafate, Sinemet SR, Elavil, Keflex, Micronase, Cardene, Fluconazole, Reglan, Wellbutrin. Allergic to deposedatives, Cipro, Reglan, Zestril.	Dopamine, Glucernia, D5, Tagamet, Dilantin, Accucheck, Procardia, Rocephin, Humulin, Nitroglycerin, Atropine sulfate, Maalox, acetaminophen, ACLS protocol, Lasix.
*HMC523	Male	52	C	Anoxic Injury	Smoker, Drinker (2-4 oz P/day)	History of high blood pressure, shunt from R-L. Intestinal problems, pulmonary fibrosis. Clubbing in extremities. Skin grafting from firev'69, Hernia surgery '85	Levophed, Dioxin, Vancomycin, Dopamine, Halcion, Ceftranxone, Lasix, Bactrum, Coumadin, Heparin, Solumdrol, Prednisone, Blood products
HMC524	Male	51	C	Subarachnoid hemorrhage	Not available. Tattoos. Positive for alcohol on admission.	Not available. Craniotomy on admission.	Heparin, Solumedrol, Dopamine, Levophed, Lasix, Mannitol, DDAVP, Ancef, Nystatin, Neomycin, Erythromycin, Phenobarbital, Nimotop, Vancomycin, Dilantin, Axid, Solucortef, Fortoz, Vitamin K, Hydrocortisone, crystalloids, colloids.

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

© DLS

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
HFC551	Female	28	C	Spontaneous cerebral bleed	Possible marijuana as teen. Smoker and drinker (1 pint Vodka/day 3-4 yrs), was in alcohol detox 4-5/95	AVM seizures '91, AVM repair. Daily meds Tegrentol since '91.	Dopamine, Mannitol, Decadron, Dilantin, Ancef, Pepcid, Pitressin, Blood Products, Crystalloids
HFC56	Female	57	C	Anoxia	Not available.	Amoxicillin, HTN, Pulmonary Fibrosis, Arthritis, Diabetes, Insulin Injections	Dopamine
*HFC586	Female	42	C	Subarachnoid Hemorrhage	Social drinker beer weekends. Moderate smoker.	Complained of hand pain while typing, tubligation 5 yrs ago. Daily meds Aspirin, T3 for sinuses	Decadron, Mannitol, Sinusitis, Digoxin, Esmolol, Pitressin, Dilantin, Veised, Ativan, Sucis, Digoxin, Dopamine, Blood Products, Crystalloids
*HMC597	Male	46	C	Subarachnoid Hemorrhage	1 pint per day social drinker x 6 yrs.	Possible exposure to Agent Orange in Vietnam. Chronic right side pain. Nov '91 fatty tumor above right knee, intertwined in muscle non-cancerous. May 1995 non-insulin dependent diabetes. Lost 30 lbs last year AMA diet. Stiff joints. Daily meds Prozac, Glucophage, Flexeril, Doxepin, Glynase.	Mannitol, Ancef, Nimodipine, Insulin
*HFC598	Female	59	C	Intracranial Hemorrhage	Non-drinker 1 PPD smoker.	History of HTN 6-7 yrs. Diabetic oral meds 1/90 insulin as of 6/90. Spasmodic stomach. History of arthritis in hands, hips, back, knees 20 yrs. Tonsillectomy 12 yrs ago. Ovarian tumor removed 1986. Orhto on knee 1988. Partial hysterectomy 19 yrs ago. Daily meds Calan, Insulin, Coumadin, Darvacet, Amytriplen, Zantac.	Ancef, Dopamine
HFC600	Female	32	C				
*HMC601	Male	41	C	Closed head injury s/p motor cycle accident	12 pack beer/week. Marijuana, cocaine (non IV) in 20s.	Laminectomy at 4 years old. Benign tumor removed 16 years ago.	Dopamine. (medical records incomplete)
HFC604	Female	42	C	Smoke asphyxiation from fire	Smoker, Non-drinker. Very active.	Bell's Palsy w/ C section pregnancy 14 yrs ago.	Dopamine, Ancef, Lasix, Calcium chloride, Potassium Bicarbonate
*HMC609	Male	48	C	Head trauma s/p Motor cycle Accident	Normal.	Positive TB test 20 years ago, treated. Broken arm as child. Allergy medicines, vitamins.	Dopamine, Epinephrine, Ancef, DDAVP, T4, KCl, insulin, Bicarb
*HFC610	Female	39	C	Intracranial Bleed	Smoker, drinker (12-18 beers/wk x 4 yrs) occasional marijuana 4 yrs ago. Military service 1974-1976.	HTN diagnosed 5 yrs ago no meds. Heart valve narrowed @ birth. Gall bladder removed 2 yrs ago. Tonsils removed.	Insulin, MgSO4, Levophed, Dobutrex, Neosynephrine, Nipride, Dopamine, Umasyn, Blood Products, Crystalloids

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

© DLS

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
*HMC618	Male	58	C	Cerebrovascular accident	Non-drinker, ex smoker	Had kidney stones 5-8 yrs ago. Broken ankle, wrist, and cracked ribs yrs ago. Ankle & appendix surgery yrs ago. Regular blood donor	Ancef, Dopamine, DDAVP
*HFC631	Female	38	C	Intracranial Bleed	Smoker, light drinker (6 pk/week)	Gallbladder removed. 3 weeks ago hospitalized with back pains	
*HFC637	Female	59	C	Cardiac Arrest	Non-smoker; Non-drinker	Saw Dr. in October for anxiety was given Psych referral. Hysterectomy 10 yrs ago. Little Arthritis. Daily meds Noratrypt	Dopamine, Dobutamine, T3 Protocol, Ancef, Blood products
*HFC674	Female	59	C	Intracranial Bleed	Non-smoker; Non-drinker	Arthritis in shoulder. Detached retina not needing surgery. Bunion removed 5 yrs ago. Bladder surgery 15 yrs ago. Daily meds Propanol, Decadron	Dopamine, NCO, Heparin, Mannitol, Sdolumedrol, Lasix, Hespal, PAU, Neasynephrine, Pavalon
*HFC685	Female	44	C	Intracranial Hemorrhage s/p fall down stairs	Secretary Signed donors card social drinker, light smoker	Good health. Daily meds Prozac.	Ancef, Solumedro, Blood Products
*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, Catapres, Tylenol w/codeine, Percocet	Dopamine, Mannitol, Decdron, Zantac, Pitressen, Succinylcholine
*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.
*HFC749	Female	74	C	CVA-ICH	Normal.	Hypertension x20 years. Congestive heart failure, pneumonia, blood infection 1 year ago. Vein surgery years ago. Arthritis. Coumadin, Vasotec, Synthroid, Retinal, Lanoxin, K-Dur, Lasix.	Heparin, Dopamine, crystalloids.
*HFC755	Female	42	C	SDH, Cerebrovascular Accident s/p aneurysm	Social alcohol use.	Hysterectomy to resolve endometriosis. Possible hypertension. Migraines x3 years. Fiorinal, Imitrex. On this admission, craniotomy.	Heparin, Mannitol, Solumedrol, Dopamine, prostaglandin, Phenergan, Compazine, Decadron, Pepcid, Cerebyx, DDAVP, Rocephan, Hespan, Pitressin, D50

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

© DLS

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
*HFC761	Female	47	C	Intracranial Hemmorage	1-2 beers/day, 10 years (heavy ETOH earlier), marijuana (quit 2 years ago). Jailed overnight, DUI, 1994.	Recovery sheet notes fibrous nodular hyperplasia in liver, else normal.	Heparin, KCl, Dopamine, Pitressin, Cardizem, nitroprusside, NTG, Ephedrine, Neosynephrine, Kefzol, Cleomycin, Norcuron, phenobarbital, colloids, crystalloids
*HMC764	Male	52	C	SIGSW	Beer and rum daily, 3-4 drinks, >20 years. Tattoos.	Hypertension, recently diagnosed, no meds. Gout in big toes. Crushed finger years ago. Laceration to leg, with large blood loss, 1989, possible blood transfusion. History of IBS. Vaccinated for Hep B.	Mannitol, Vecuronium, Centanyl, Versed, tetanus, T4, Solumedrol, Dopamine, D5, crystalloids, Hespan, Heparin, Dobutamine, Ancef, Lasix
*HMC782	Male	57	C	Cyanide OD, cardiac arrest	Social alcohol	Family history of cardiac disease; cardiography notes CAD. Kidney stones x3 years. Lyme vaccine (2 of 3 shots). High cholesterol. History of hives, shingles. Eye infections 6 years ago from contacts. Lipitor, Proscar, Flonase, vitamins, fish oil.	Charcoal, sodium nitrate, sodium thiosulfate, Levophed, Vasopressin, Labetalol, Kphos, T4, D5W.
*HFC783	Female	74	C	CVA	Beer a month.	Hypertension, on meds for 3 years. Broke ankle 15 years ago. Lived near uranium mine. Craniotomy on admission.	Mannitol, Dopamine, Digoxin, Dilantin, Cardizem, Phenigran, Labetolol, D5, blood products, albumin.
HFC791	Female	67	C	Stroke	Smoker, Beer on holidays only.	Burned back '56. Goiter removed '55. Motor vehicle accident '55. Broke leg 30 yrs ago. Allergic to penicillin.	DDAVP, Solumedrol, Ancef, Dopamine
HFC794	Female	47	C	Cerebrovascular accident	Smoker, rare social drinker	Partial hysterectomy for cancer in early 1980's. Frequent constipation. Left broken toe. Edema as a child	
*HFC803	Female	21	C	Head trauma s/p Motor Vehicle accident	Non-smoker, Alcohol with meals on trip to Europe 2000. Pierced ears and naval both professionally done.	Broken collar bone and pelvis at age 3-4. Jaundice at birth.	Dopamine, T4, TPN
*HFC815	Female	49	C	Head trauma, TBI s/p MCA	Smoker, light drinker one beer a day.	Famvir (1 BID). Tubal libation 1989. WBC 11.19, temp 95-102, probably neurogenic. On admission 6/15/03, SGOT 38 ALT 25. Final on 6/16/03 SGOT 1305, ALT 729, T Bili 1.1, Direct 0.4. Liver enzymes unusually high.	Dopamine, Epinephrine, Neosynephrine, Dilantin, Rocephin, Protonix
*HFC817	Female	56	C	Subarachnoid Hemmorage	Non-smoker; Non-drinker	Broken left wrist '93. High blood pressure since '92 Arthritis in right knee. Daily meds Dioran, Valium	Pe4pcid, Labetolol, Insulin, Ancef, "Nimodipine, Dulcolax strup, Neutra, Dilantin, Tylenol syrup, MS,
HFC820	Female	62	C	CVA-ICH	3 glasses wine daily	Hiatal hernia; biopsy revealed fibrous tumor. Hypertensive x3 years, treated with meds (Diovan, Tricor). Cellulitis right leg summer 2003. Colon polyps removed x3-4 years (2-3 nonmalignant polyps, repeat colonoscopy clear). Allergic to sulfa drugs.	Mannitol, KCl, Solumedrol, Dopamine, Labetolol, morphine, Tylenol, Diprivan, Norcuron, Cefazolin, Versed, Pepcid, crystalloids

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

© DLS

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
*HFC823	Female	48	C	Anoxia s/p MVA	1 glass wine every 6 months	Vitamins, Prozac. Received Hep B vaccination. 2 episodes of down time, x4 min and x3 min. Lacerations on liver from MVA.	T4, Dopamine, Epinephrine, Levophed, Ancef, Diflucan, Versed, Pavulon, Heparin, Mannitol
HFC828	Female	28	C	Anoxia s/p CVA	2-4 beers a week. Jail 10 years ago, overnight.	Asthma x2 years, treated with Albuterol. MS x2 years, treated with Copaxone. Kidney stones. 6 months post-partum. Exposure to chemical spill in photo developing lab 3 years ago. Broken arm at 5 years old. Screening record notes left lobe normal, right lobe central lobular necrosis.	Heparin, Dopamine, albumin, Regitine
*HFC830	Female	32	C	Intracranial Hemorrhage	Smoker, light drinker 1 drink month or less.	History of seizure, last one 11 years ago. Aneurysm at 9, surgery at 14, C section 1993. Kidney and yeast infection treated with ATB. UTI x1 yr ago. Neurontin, Verapamil (Verapamil?).	Heparin, KCl, Levophed, Vasopressin, Phenylephrine, Fentanyl, blood products, albumin, NS
*HFC833	Female	52	C	Cerebrovascular accident	Drank once every 10 days x5 years. No IV drugs; used cocaine, smoked marijuana once a week, drug history over 5 years.	Hysterectomy 5 years ago, acid reflux x2 years, hypertension x2-3 years. Albuterol, Prilosec, Lotrel.	KCl, Dopamine, Vasopressin, Labetolol, Protonix
*HFC835	Female	52	C	Cerebrovascular accident	Non-smoker; Non-drinker	No past medical history. Had cold with sinus infection, temporarily using steroid inhaler.	Heparin, T4, Dopamine, Levophed, albumin
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, Cefepine, Solumedrol, Humulin, Topramycin, Flagyl. (records unclear)
*HMC856	Male	51	C	Head trauma s/p fall.	Scotch or wine 4-6 drinks a day. Marijuana use in his 20s (x30 years ago).	Aortic stenosis at birth; 2 surgical repairs at age 10 and 28. Possible hypertension per NOK. Hematomas on legs, possibly due to Coumadin per PCP. Probably tonsillectomy. Periodic migraines. Post fall, refused to seek treatment, took 2 hydrocone, bottle of wine and 2 gins night before admission to hospital.	Narcan, Adenosine, Versed, Neosynephrine, Levophed, Vasopressor, Dopamine, Cardizipem, KCl, T4, Pitressin, Insulin, Ancef, Solumedrol, Accucheck, Labetalol, Kphos, MgSulfate.
HFC863	Female	46	C	Anoxia	Drug use (crack cocaine, marijuana). Jail time - 6 months, 1998, drug charges.	Asthma, depression. Flovent.	Heparin, Levophed, T4, insulin, crystalloids, albumin.
HFC865	Female	53	NA	Head trauma s/p MVA.	Occasional wine.	Kidney stones a few years ago. Possible liver hemangioma.	Heparin, Regitine, Versed, blood products

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

© DLS

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
*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.
*HMC906	Male	67	C	ICH-Stroke	Alcohol; 2 drinks or less per month.	Normal healthy	Dopamine, Nicardipine, Levothyroxine, NS, Insulin, Lasix, Zosyn, Vancomycin, Labetolol, Hydralazine, Fentanyl, Mannitol, Crystalloids, Heparin, and Pancuronium. 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.
*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.
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.
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.

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

© DLS

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
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
*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.

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

© DLS

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
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, Neosyneprine, 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.
*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, Zoloft, Ibuprofen, Nerve medicine for back (as per NOK).	Epinephrine, Narcan, KPhos, Labetalol, KCl, MgSO ₄ , Fluconazole, Ancef, Vasopressin, Levophed, Neosyneprine, 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.

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

© DLS

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
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.
*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.

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

© DLS

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
*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.
*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.

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

© DLS

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
*HFC954	Female	32	C	Head Trauma 2 nd Blunt Injury	ETOH (Beer, Wine) 2 x / wk X 10 yr.	No meds taken at home (as per NOK).	Levophed, Dopamine, Neosynephrine, Solumedrol, Lasix, T4, Vasopressin, Narcan, Pepcid, Combivent, Zosyn, Vancomycin, CACI, KCI, Betadine, Amphotericin, Epinephrine, Norepinephrine, Heparin, Fentanyl, Propofol. 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.
*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.

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

© DLS

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
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.
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.

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

© DLS

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
*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.
*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
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.
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, Solucortef, Cefazolin, MgSulfate, Propranolol, DDAVP, Dilantin, D5W, Pepcid, Vecuronium, Midazolam, Fentanyl, Succinylcholine, Etomidate, Senna, Heparin. Flushed and stored in SPS-1.

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

© DLS

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.
*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.
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, NaHCO3, Propofol, Levophed, Insulin, Albuterol, NS, Solumedrol, Kayexalate, Hespan, CaCl, Heparin, Mannitol, Lasix. Flushed and stored in UW solution.
*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.

© DLS

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.
HMC1030	Male	32	C	Anoxia			Dopamine, T4, Insulin, Flushed and stored in HTK.
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.
*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, Kepra, Sodium BiCarbonate, LR, Ancef, KCl, KPhos, Flushed and stored in SPS-1.
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.

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

© DLS

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
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.
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.
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, Zolof until June 20115, MVI, Glucosamine (as per NOK).	Levophed, Albuterol, DuoNeb, Precedex, Prednisone, Solumedrol. Lasix, Bumex. Flushed and stored in HTK.

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

© DLS

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
*HMC1071	Male	50	C	Anoxia 2nd 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.
*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, HLD. 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,

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

© DLS

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.
- Treat all products containing human-derived materials as potentially infectious, as no known test methods can offer assurance that products derived from human tissues will not transmit infectious agents. 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.



Quality Assurance

10 September 2024

Date

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

© DLS