

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

# UltraPool™ HLM 150

Catalog No.	452117
Lot No.	38300
Qty./Package	0.5 mL
Protein Content	20 mg/mL in 250 mM sucrose
Storage Conditions	Store at -80°C
Release Date	2024 April
Expiration Date	2033 November

## Assay Results

Enzyme Measured	Assay	Enzyme Activity [in pmol/(mg x min)]
Total P450	Omura and Sato	320 pmol/mg
OR	Cytochrome c Reductase	150 nmol/(mg x min)
Cytochrome b <sub>5</sub>	Spectrophotometric	380 pmol/mg
CYP1A2	Phenacetin O-deethylase	420
CYP2A6	Coumarin 7-hydroxylase	770
CYP2B6	Bupropion hydroxylase	260
CYP2C8	Paclitaxel 6 $\alpha$ -hydroxylase	140
CYP2C9	Diclofenac 4'-hydroxylase	2700
CYP2C19	(S)-Mephenytoin 4'-hydroxylase	51
CYP2D6	Bufuralol 1'-hydroxylase (The amount of activity inhibited by 1 $\mu$ M quinidine.)	93
CYP2E1	Chlorzoxazone 6-hydroxylase	2200
CYP3A4	Testosterone 6 $\beta$ -hydroxylase	4100
CYP4A11	Lauric acid 12-hydroxylase	1700
FMO	Methyl p-Tolyl Sulfide Oxidase	3700
UGT1A1	Estradiol 3-Glucuronidation	1700
UGT1A4	Trifluoperazine N- Glucuronidation	690
UGT1A6	Serotonin Glucuronidation	10000
UGT1A9	Propofol Glucuronidation	2800
UGT2B7	AZT Glucuronidation	1400
<b>Western Blot Results</b>		
CYP3A4	Western Blot	92 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 MgCl<sub>2</sub>, 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 MgCl<sub>2</sub>, and 25  $\mu$ 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](mailto: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
HFC25	Female	66	C	ICH	Tobacco use.	HTN, Stomach ulcer, Rheumatoid Arthritis, Immodium, Zinc Teap, Augmentin, Diflucan, Vancuron, Celebrex, Nevantin, Furosemide	Not available.
*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, Levothyroxine, Levo drip and Narcan
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
*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.
*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.
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.
*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.

**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](mailto: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
*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. Zydys, Lovastatin, Acetaminophen, Aspirin, Bupropion, Dicyclomine, Escitalopram, Flurazepam, Gabapentin, Lisinopril, Metformin, Ziprasidone (as per NOK).	NS, Dopamine, Levaquin, Clindamycin, Insulin, Levofloxacin, Narcan, Norcuron, Pitressin, Potassium Chloride, Sodium Bicarbonate, Solumedrol, Water, Albuterol, Atrovent, Magnesium.
*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, Levofloxacin, Vasopressin, Albuterol, Levophed, Insulin, Dopamine, 0.45 NS, Levofloxacin, D5W, Vasopressin, PRBC, 5% Albumin, KCl in NS, D5, Mannitol, T4, and Crystalloid. Flushed and stored in UW solution
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.
HFC477	Female	52	C	ICH-Stroke	Cigarettes: 1/2 PPD X 2 months. 2 PPD X past 30-40 yrs. Beer: Up to 10 beers, 4-5 times a week X 20+ yrs. 10 beers a day only X past 10 years.	Stroke. Diagnosed with Hypertension 20+ yrs prior and treated with Lisinopril. Seizures. Diverticulitis and Ulcers. Kepra, Lortab, Xanax, Lisinopril, and Aspirin (as per NOK)	Nicardipine, Mannitol, Solumedrol, Albumin, 0.45 NS, Dopamine, Ancef, PRBCs, Nipride, Crystalloids, Heparin, Thorazine, Fentanyl, Rocuronium and Vasodilators. Flushed and stored in HTK.
*HFC499	Female	28	C	Head Trauma-2nd 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,
*HFC522	Female	58	C	Not available	Smoker, light drinker 1-2 drinks week.	Treated for Asthma. Spoken with doctor concerning the possibility of heart disease. Suffered from arthritis in both hands.	Procardin, Ecotrin
*HFC523	Female	51	C	Anoxia-2nd 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.
*HFC586	Female	42	C	Subarachnoid Hemorrhage	Social drinker beer weekends. Moderate smoker.	Complained of hand pain while typing, tubigation 5 yrs ago. Daily meds Aspirin, T3 for sinus	Decadron, Mannitol, Sinusitis, Digoxin, Esmolol, Pitressin, Dilantin, Veised, Ativan, Sucis, Digoxin, Dopamine, Blood Products, Crystalloids
HFC589	Female	48	C	Aneurysm	3-4 whiskeys per day x10 years. Tobacco	Tubal ligation 18 years ago. Liver inflammation. Menopause, night sweats x1 year. IIM screening form notes findings consistent with chronic hepatitis, possibly from alcohol, autoimmune or viral (note: recovery sheet notes liver looks fine, serologies and PCR results are negative.) Possible Hep B vaccination while EMT. Antibiotics 6 weeks ago, abscessed tooth. Allergic to penicillin.	Heparin, Mannitol, Lasix, Solumedrol, Dopamine, Neosynephrine, Pitressin, Nipride, Dilantin, T4, Vascomycin.
*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				
*HFC613	Female	33	C	Anoxia s/p seizure	Social alcohol	Diabetes, noncompliant with insulin x6 months, Glyburide sporadically. Tendinitis. Appendectomy 1977. Obesity. Hypertension. Chronic poor health. Ovary infarcted (tube twisted). Digestive problems associated with headaches. Otitis inflammation. ASA. Reflex for recent ear infection. Allergic to strawberries.	Heparin, Mannitol, Dopamine, Dobutamine, Narcan, Dilantin, Cigforan, Ampicillin/Sulbactrim, Verapamil, Lasix, albumin, crystalloids, colloids.
*HFC631	Female	38	C	Intracranial Bleed	Smoker, light drinker (6 pk/week)	Gallbladder removed. 3 weeks ago hospitalized with back pains	
*HFC640	Female	67	C	Cerebrovascular intracranial hemorrhage	Smoked <pk day quit 20 yrs ago. 1 glass red wine at night.	HTN 10 yrs on medicine (Catapril) daily Zantac. Hysterectomy 20+yrs ago. Flu shot 12/96.	Dopamine, Pitressin, Mgs04, Heparin, NAHCO3, Pavulon, Cace, D50, Fentanyl, 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](mailto: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
*HFC650	Female	58	C	Cerebrovascular Accident, Intracranial Bleed s/p Aneurysm	Social alcohol use.	Hypertension. High cholesterol. Benign bilat breast biopsies. Peptic ulcer late 70s. Arthritis in feet, knee stiffness s/p MVA 1983. Beroca Plus, Tylenol with codeine, Zantac, Tenoretic, Vasotec, Zocor. Asbestos exposure 10 years ago.	Heparin, Solumedrol, Regitine, Neosynephrine, Dopamine, Mannitol, Ancef, Zantac, MgSO4, CaGluconate, Dilantin, albumin, crystalloids.
HFC665	Female	64	C	Cerebrovascular Accident, Intracranial hemorrhage	Normal.	Early stage emphysema, 5/96. CHF from pneumonia, 5/96, treated with Lanoxin for atrial fib. Diabetes x10 years, controlled with Glucotrol (oral); 5/96 became IDDM. Hypertension x19 years. Diverticulitis. Hysterectomy, 1965. Cholecystectomy, 1976. Lasix, Tegretol, MicroK, Chlorprompemde.	Heparin, Dopamine, D5, Nipride, Ancef, Lasix, crystalloids.
*HFC666	Female	52	C	Intracranial Hemorrhage	Social alcohol use.	Cholecystectomy 1972. Immune Thrombocytopenic Purpura diagnosed 1993, in remission since 1994. Angina recently diagnosed. Esophageal reflux Nitroglycerine, Prilosec, Premarin.	Heparin, Fentanyl, morphine, Demerol, Valium, Versed, Pitressin, Nimodipine, Dopamine, insulin, Decadron, Cimetidine, Phenergan, albumin, D5, Ancef, Mg, Phos, Calcium, Lidocaine, Tagamet, DDAVP.
*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
*HFC677	Female	63	C	Subarachnoid hemorrhage	Social alcohol use.	Borderline diabetic x6 years. Coronary bypass 1994. Stent 2 years ago. Hypertension >10 years. Arthritis in hands and hips. Hysterectomy, cholecystectomy, cataract removal with implants. Coumadin, Atenolol, Imdur, Axid, Zestoretic.	Heparin, Zemuron, Dopamine, Dobutamine, Pitressin, crystalloids, colloids, Cerebyx, T3, insulin, Solucortef, D50W, NaBicarb, MgSO4, CaChloride.
*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
*HFC696	Female	45	C	Closed head injury s/p motor vehicle accident	Normal.	Jaundice 15 years ago. Malaria 15 years ago, contracted s/p transfusion while treating jaundice; recurrence 7 years ago. Hysterectomy 10 years ago. Tetanus shots 16 years ago to treat animal bites. Kidney stones 1 year ago. Benign skin lesions removed from face >1 year ago. Lived in Africa for 15 years, last there 6 years ago.	Dopamine, Pitressin, DDAVP, Lidocaine, Folic acid, Selenium, Cenolate, Solumedrol, Mannitol, Heparin, CaGI, blood products
HFC703	Female	29	C	Overdose	Social alcohol. Cocaine (inhaled), crack x5 months 2 years ago. Positive for amphetamines, cocaine on admission, drug history denied by family, confirmed by boyfriend. Possible track marks.	Possible abortions, 3/97, 11/97. Possible recent infection with chlamydia. Oral contraceptives.	Heparin, Dopamine, crystalloids, Cardizem, Solucortef, NAHCO3, D5, Ticarcillin, albumin, blood products, Epinephrine, Narcan
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
*HFC817	Female	56	C	Subarachnoid Hemorrhage	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, Peppid, crystalloids
*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

**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](mailto: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
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)
*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, Atacurium, Mannitol, Labetalol, Pepcid, Ancef, Dopamine, Remifentaniil
*HFC903	Female	60	C	Anoxia-2nd 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 Lvxyyl.(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 2nd 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.
*HFC909	Female	64	C	Anoxia 2nd 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.
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.
*HFC931	Female	19	C	Anoxia 2nd 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.
HFC934	Female	54	C	CVA 2nd 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, Zvox, Cefepime, Levothyroxine, Lasix, KCL, Furosemide, Zithromax, Rocephin, Propofol, Heparin, Magnesium Sulfate. Flushed and stored in HTK.
*HFC944	Female	44	C	CVA 2nd 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 2nd 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](mailto: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
HFC952	Female	40	C	Anoxia 2nd 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.
*HFC954	Female	32	C	Head Trauma 2nd 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, KCl, Betadine, Amphotericin, Epinephrine, Norepinephrine, Heparin, Fentanyl, Propofol. Flushed and stored in UW solution.
HFC955	Female	66	C	CVA 2nd 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, Metoprolol, Cardene, and Vasopressin. Flushed and stored in UW solution.
HFC957	Female	36	C	CVA 2ND 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.
*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, Impipenem, NS, Diuretics, T4, Vasodilators and Antihypertensives. Flushed and stored in SPS-1.
HFC963	Female	74	C	CVA 2ND 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 2nd 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 2nd 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.
*HFC966	Female	65	C	CVA 2nd 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 2nd 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.

**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](mailto: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
*HFC970	Female	75	C	CVA 2nd 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.
*HFC979	Female	27	C	Head Trauma 2nd 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.
*HFC982	Female	62	C	CVA 2nd 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.
*HFC987	Female	63	C	CVA 2nd 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)	Labeltolol, Colace, MgHydroxide, Dulcolax, Novolog, D50W, Artificial Tears, Protonix, Combivent, Cardene, KCl, Zofran. Flushed and stored in UW solution.
*HFC993	Female	65	C	Anoxia 2nd Cardiovascular	Cigarettes- 2 PPD x 45 yrs. ETOH (Beer, Liquor) a 5th 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.
*HFC997	Female	48	C	CVA 2nd 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 PMH: HLD, hypothyroidism.	Dopamine, Levophed, Famotidine, Kepra, Vasopressin, Dextrose, T4, Insulin, DDAVP, Heparin. Flushed and stored in HTK.
HFC1005	Female	49	C	Anoxia 2nd 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.
HFC1009	Female	24	C	HT 2nd 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, Kepra, Mannitol, KPhos, NaPhos, Heparin. Flushed and stored in HTK.
*HFC1013	Female	40	C	CVA 2nd 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.

**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](mailto: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
*HFC1017	Female	63	C	Other 2nd 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.
*HFC1021	Female	46	C	Pneumonia 2nd 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.
*HFC1028	Female	46	C	CVA 2nd 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.
*HFC1035	Female	59	C	CVA 2nd 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, Lisinpril, Pregabalin, Bupropin (meds taken at home as per NOK)	Esmolol, Solumedrol, Insulin, Antihypertensives, DDAVP, KCl, Labetalol, Nipride. Flushed and stored in SPS-1.
*HFC1043	Female	53	C	Anoxia 2nd 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 2nd 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
*HFC1058	Female	47	C	Anoxia 2nd 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, Lantus, Januvia.	Levophed, Vasopressin, Insulin, Propofol, Zosyn, Decadron, Heparin. Flushed and stored in SPS-1.
*HFC1067	Female	56	C	Anoxia 2nd 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.
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 Ablation 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, Neosynephrinemm and Solucortef. Narcan and Romazicon administered
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.

**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](mailto: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
HMC274	Male	37	C	Anoxia	Cigarettes; 1 1/2 PPD x 15 yrs. Wine; once in a while; not a heavy drinker. Heroin; snorted on and off for years; last used 2 weeks ago. Upon admission pt was positive for opiates and Benzos.	Pt had pacemaker inserted in 20002, Heartburn; lost 15 lbs over last 2 wks d/t detox. Multivitamin, Beritol, B12, Trazadone, Lasix, CaCl, NS and Levophed.	Epi and Atropine.
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.
HMC278	Male	26	C	Cerebral Edema	Cigars; Every other day or less x 1 yr. Beer and Whiskey; Socially since college.	Hypertension; Diagnosed 12 yrs ago. Diet controlled. Multivitamins and Metamucil (as per NOK).	Dopamine, Neosynephrine, Levothyroxine, Vasopressin, Ancef, Fortaz, Crystalloids, Coloids, 5% Albumin, Heparin, Mannitol, T4 and Insulin. Flushed and stored in Celsior.
HMC282	Male	41	C	ICH-Stroke, MVA	Tobacco- 1/2 PPD X 25 yrs. Beer- 1/ mo since age 21. Cocaine/Heroine- 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.
*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, Zoloft, and CHF meds (as per NOK).	Acetaminophen, Acetylcysteine, Calcium Gluconate, Dextrose, digoxin, Epinephrine, Esmolol, Furosemide, Heparin, Hydralazine, Levetricetam, Mannitol, Metoprolol, Neosynephrine, Piperacillin, Vancomycin, NS, Levophed, Solu-Medrol, Zosyn, Albuterol, Insulin, Lasix and Vecuronium. Flushed and stored in Custodial.
*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.
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, Cystalloids, Levophed, and Heparin. Epinephrine and atropine were administered. Flushed and stored in UW solution.
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.
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.
*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.
*HMC509	Male	44	C	Closed head injury motorcycle vs car.	Smoker, drinker (a lot mostly beer x 25 yrs).	HTN borderline never prescribed meds.	Dopamine, lidocaine, Epinephrine, Atropine, Dilantin, Zantac, Rocephin, Kefzol, Calcium chloride, MS04, DDVAP, KC1
*HMC510	Male	22	C	Head Trauma-2nd 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.
*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

**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](mailto: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
HMC542	Male	32	C	Closed head injury s/p motor vehicle accident	Light smoker social drinker	Healthy gave birth 2 yrs ago	Dopamine, Ancef, Albumin, Blood Products, Crystalloids
*HMC552	Male	47	C	Brainstem Stroke	Smoker, quit drinking 14 yrs ago. Marijuana 1 x 2-3 yrs ago	Cramps in legs. Daily meds Tegrentol	Dopamine, Pitressin, Ancef, Heparin, K Phos, Carafate, Zantac, Solumedrol
*HMC562	Male	54	C	Cerebrovascular accident	Normal.	Hypertension. CAD, 3 previous MIs. Limited patient history available.	Dopamine, Dobutamine, Dobutex, Hespan, Levophed, Lidocaine, NaPhos.
*HMC568	Male	32	C	Anoxia, CPA	6 pack beer/day. Occasional marijuana.	Hypertension, chest pain x10-15 years, complaint with meds. Tenonmin, occasional Valium.	Inocor, Levophed, Dopamine, Dobutamine, Tenonmin, occasional Valium.
HMC571	Male	44	C	Subarachnoid hemorrhage	Up to 12 beers/day. Marijuana. Possible IV drug use. Jail, 60 days, 5 months ago for disorderly conduct. Pierced ear, nonprofessional.	Possible tonsillectomy. Right leg injury 1972. Hematuria 2 weeks ago.	Heparin, Diazepam, Pavulon, DDAVP, Digoxin, Morphine, Thiamine, Valium, Zantac, Decadron, Mannitol, Dilantin, DTPA, Dopamine, Dobutamine, Lidocaine, Nipride, D5, albumin, crystalloids.
*HMC575	Male	66	C	Cardiac arrest	Occasional alcohol; heavy alcohol use >5 years ago. Tobacco	Arthritis in back and hips.	Heparin, Dopamine, Dobutamine, Lasix, Lidocaine, Riopan, insulin, Ancef, Zantac, bicarb, albumin, crystalloids.
*HMC635	Male	66	C	Intracranial Hemorrhage	Social alcohol use.	Cardiomyopathy. Hypertension x3 years. Appendectomy at 9 years old. Treated 4-5 years ago for irregular heart beat. Chronic renal insufficiency. Micro K, Coumadin, Lasix, Lanoxin, Cipro, Procardia.	Heparin, Pavulon, Dopamine, crystalloids.
HMC668	Male	39	C	Intracranial Hemorrhage	Social alcohol, none in past year. Occasional marijuana. Tattoos, professional.	C section 9 years ago. Right elbow surgery 3-4 years ago. Arthroscopic surgery, both knees, 12-13 years ago. Frequent bronchial infections. Hypertension x3 months, treated with Procardia. Aneurysm clipped after admission.	Heparin, Mannitol, Dopamine, Nipride, Ancef, Thyroxine.
HMC724	Male	49	C	Cerebrovascular Accident s/p aortic dissection	Rarely drank. Tattoo 2 years ago. Jail, 1 night, when 16.	AAA repair, valve replacement, on admission 6/24/99. Meningitis, 1965.	Dilantin, Lidocaine, Versed, succinyl choline, Pavulon, Mannitol, Cefazolin, Neosynephrine, insulin, Decadron, blood products, crystalloids, Dopamine.
HMC812	Male	22	C	Anoxia s/p hanging suicide	Smoker, Heavy drinker at least a 6 pk /per day. Used cocaine and marijuana. No IV drug usage. Had tattoos not professionally done spent 3 days jail in 2001.	Healthy	Dopamine, Neopride, Insulin, Mannitol, KCl, MgSO4, Abcef, NaHCO2, Tylenol, NEO, EPI, Crystalloids
*HMC911	Male	53	C	CVA 2nd 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.
*HMC913	Male	37	C	Anoxia 2nd 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.
HMC922	Male	42	C	Other 2nd 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, Lovexox, Ativan. Flushed in Custodial. Stored in HTK.
HMC928	Male	37	C	ANOXIA 2nd 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.

**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](mailto: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
HMC930	Male	56	C	Anoxia 2nd Electrocuted	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
*HMC933	Male	39	C	Head Trauma 2nd 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.
*HMC936	Male	60	C	Head Trauma 2nd 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, Neosynephrine, 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 2nd 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, MgSO4, Fluconazole, Ancef, Vasopressin, Levophed, Neosynephrine, Levothyroxine, Solumedrol, Mannitol, Heparin. Flushed and stored in SPS-1
HMC938	Male	54	C	Anoxia 2nd 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 2nd 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 2nd 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 2nd 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.
*HMC947	Male	68	C	CVA 2nd 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.

**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](mailto: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
HMC949	Male	50	C	Anoxia 2nd 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.
*HMC950	Male	50	C	Head Trauma 2nd 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.
HMC953	Male	76	C	CVA 2nd 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.
*HMC960	Male	57	C	ICH 2nd 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.
*HMC962	Male	60	C	CVA 2nd 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, Solucortef, T4, Insulin, Heparin, Mannitol , Rocuronium, Dexamethasone. Flushed and stored in SPS-1.
HMC965	Male	46	C	Anoxia 2nd 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.
HMC974	Male	33	C	HT 2nd 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 2nd 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 2ND 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, Numbutal, Magnesium Sulfate, Cefepime, Vancomycin, Docusate, Famotidine, Versed, Etomidate, Vecuronium and Platelets. Flushed and stored in SPS-1
HMC988	Male	52	C	Anoxia 2nd 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.

**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](mailto: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
HMC992	Male	30	C	HT 2nd 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 2nd 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.
*HMC1003	Male	32	C	Other 2nd 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 2nd Drug Intoxication	Cigarettes: 1.5 PPD since 18 yrs old. ETOH: (Distilled Spirits) Drank unknown amount for unknown duration. Heroine: 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.
*HMC1007	Male	61	C	HT 2nd 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 2nd 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, Hesperin, 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. Liposil, 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.
HMC1027	Male	66	C	Head Trauma 2nd 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, Benztrapine 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.

**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](mailto: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
HMC1030	Male	32	C	Anoxia			Dopamine, T4, Insulin, Flushed and stored in HTK.
HMC1036	Male	56	C	HT 2nd 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 2nd 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.
HMC1048	Male	36	C	CVA 2nd 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 slowing 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 2nd 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 2nd 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 2nd 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 2nd 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 2nd 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.
*HMC1064	Male	66	C	CVA: 2nd to natural causes	Tobacco use: 1/2 ppd x 20 yrs, quit 10 yrs prior to death	Hypertension: diagnosed 10 yrs prior, compliant w/ meds, Diabetes: diagnosed 10 yrs prior, compliant w/meds, GI disease: IBS over last 2 years - no treatment, colonoscopy showed few polyps( benign), MVI, Methocarbamol, Coumadin, Lisinopril, Hydrochlorothiazide, Atenolol, Sotalol, Metformin, Tramadol, Rosuvastatin	Vecuronium, Heparin, Levophed, Solumedrol, Lasix, T4, Insulin, Vasopressin, Ancef, Albumin
HMC1066	Male	46	C	Anoxia 2nd 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.

**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](mailto: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
HMC1070	Male	55	C	Anoxia 2nd 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 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](mailto: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.



08 May 2024

---

**Quality Assurance**

**Date**

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

© DLS