

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

# Gentest<sup>®</sup> UltraPool<sup>™</sup> Human S9 150

Catalog No.	452116
Lot No.	2403281
Qty./Package	1.0 mL
Protein Content	20 mg/mL in 150 mM KCl, 50 mM Tris (pH 7.5), 2 mM EDTA
Storage Conditions	Store at -80°C
Release Date	2024 September
Expiration Date	2034 April

## Assay Results

Enzyme Measured	Assay	Enzyme Activity [in pmol/(mg x min)]
CYP1A2	Phenacetin O-deethylase	130
CYP2A6	Coumarin 7-hydroxylase	260
CYP2B6	Bupropion hydroxylase	58
CYP2C8	Paclitaxel 6 $\alpha$ -hydroxylase	61
CYP2C9	Diclofenac 4'-hydroxylase	590
CYP2C19	(S)-Mephenytoin 4'-hydroxylase	7.3
CYP2D6	Bufuralol 1'-hydroxylase (the amount activity inhibited by 1 $\mu$ M quinidine)	19
CYP2E1	Chlorzoxazone 6-hydroxylase	600
CYP3A	Testosterone 6 $\beta$ -hydroxylase	960

- All cytochrome P450 assays were conducted at 1.6 mg/mL protein (except CYP2B6 which was at 0.2 mg/mL, CYP2C19 which was at 0.6 mg/mL, and CYP3A which was at 1 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 and CYP3A 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 pH 7.5 was used for CYP2A6 and CYP2C9). Activities expressed as pmole product per (mg protein x minute).
- 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
- The pool is comprised of equal milligrams of S9 per donor.
- HAZARD WARNING:** This S9 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.
- Donors with CMV serology unknown are identified with a double asterisk. Donors CMV negative for serology are unmarked.
- Safety Recommendations:** Safety assessment indicates this product is hazardous; refer to SDS [Safety Data Sheet] for safety information. Handle in accordance with good industrial hygiene and laboratory safety practices.

*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/>



18 September 2024

---

**Quality Assurance**

---

**Date**

**Donor Information**

Specimen	Gender	Age (years)	Race	Cause of Death	Social History	Medical History	Medication given during Hospitalization
*HMC178	Male	41	C	CVA-Stroke	Meth, wine and occasional vodka, 4 x week, 1-2 glasses x 5-6 yrs.	Crohn's Disease, cystic acne on face, Prozac.	Dopamine, Neosynephrine, Thyroxine, Vasopressin
*HMC225	Male	21	C	Head Trauma-MVA	Cigarettes; 1 PPD x 5 yrs. Occasional binge alcohol since age 9. Marijuana; stopped in 2004. Cocaine and mushrooms stopped in 2003.	healthy	Vancomycin, Levaquin, Vasopressin, Insulin, Levophed, Ancef, Cefamine, and TPN.
HFC227	Female	46	C	CVA-Stroke	Non-smoker; Non-drinker	Asthma: Diagnosed at age 5. Enbrel, Folic Acid, Iron and Nipride (As reported by NOK) Patient had plastic surgery; breast augmentation, liposuction on thighs/abdomen/upper arms and tummy tuck. Head CT showed subdural hemorrhage, which was possibly old, and pt had 2 month history of severe headaches.	Levophed, Dilantin, Dopamine, Epinephrine, Dobutamine and T4. Flush and Storage in UW solution.
HMC230	Male	21	C	Head Trauma	Beer and Liquor; unknown quantity on weekends x 2-3 yrs.	healthy	Heparin, Refitine, Vasodilators, and Morphine.
*HFC233	Female	37	C	Anoxia 2 <sup>nd</sup> Cardio respiratory	Cigarettes- 1 1/2 PPD past 10 years. Quit 1-2 yrs ago. Mixed drinks- 1 q wk. Drug use in adolescence but none for last 25-20 yrs. Drug screen was positive for Methadone, Marijuana, and Propoxyphene.	Hypertension- Diagnosed 5 yrs prior. Klonopin, Diazepam, Vicodin, Neurontin, Soma, Lisinopril, Estratab, and Ultram (as per NOK)	Dopamine, Lopresser, Clonidine, Vasolec, Solumedrol, Vasopressin, Rocephin, and Celepime. Flushed with UW solution.
*HFC234	Female	23	C	Anoxia 2 <sup>nd</sup> suicide	Non-smoker; Non-drinker		Versed, Fentanyl, Levophed
*HFC241	Female	38	C	CVA	Non-smoker; Non-drinker	Suffered Migraines and tremors of her hands and face. GERD. Multivitamin, Levaquin, Propranolol, Nortriptyline, Folic, Methotrexate, Prevacid, meloxicam, Bultalbital/APAP/Caffeine and Metoclopramide (as per NOK).	Multivitamin, Levaquin, Vacomycin, T4, Zosyn, Maxipime, Cleocine, Cancidas, Primaxin, Magsulfate, Acyclovir, Calcium Gluconate, Prtotonix, Ampicillin, KCL, Propanol, Nortriptyline, Heparin, Bactrim, Solumendrol, Diflucan, Hydrocortisone, FREE H2O, Vasopressin, Dobutamine, Lasix, Folic, Methotrexate, Prevacid, Meloxicam, Ativan, Butalbital and Metoclopramine.

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

© DLS

**Donor Information**

Specimen	Gender	Age (years)	Race	Cause of Death	Social History	Medical History	Medication given during Hospitalization
HM242	Male	58	AA				
*HMC243	Male	52	C	Head Trauma	Chewing Tobacco X 38 yrs. Hard liquor X 35 yrs; Beer- more recently; donor was alcoholic.	Pneumonia diagnosed 2-3 months prior. Hypertension; diagnosed 20 yrs ago; compliant with meds. Acid Reflux. Kidney stones; 8-10 yrs ago and 20 yrs ago. Crestor and Verapamil (as per NOK).	Levophed, Dopamine, and Propofol. Flushed and stored in HTK.
HFC248	Female	71	C	Head Trauma	Cigarettes- 1 PPD x 20 yrs. Wine- 3 glasses daily. Vodka-occasionally. Upon admission ETOH was 188.	Mini-stroke one year prior. Treated for HTN for 10-12 yrs; compliant with meds. Seizure; compliant with Keppra. IDDM since age of 30; compliant with Lantus injections. Diagnosed 12 yrs ago with IBS; compliant with Prilosec. Stent place in kidney 5 yrs prior. Keppra, Lantus injection and Prilosec (as per NOK).	Neosynephrine, Levohed, Pepcid, Senokot, Kondrumel, Vitamin K, NaHCO3, Hydrocortisone, Insulin, and T4. Flushed and stored in HTK.
*HMC249	Male	60	C	CVA	Cigarettes; 3-4 PPD x 40 yrs. Beer; 3 x 3 night q week.	Not available.	Nimodipine, Protonix, Dopamine, Vasopressin, Amiodarone, Hespain, Neosynephrine, Mannitol, Lasix, Heparin, Thorazine, and Insulin. Flushed and stored in UW soln.
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, Neosynephrinem and Solucortef. Narcan and Romazicon administered
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, Amoxicillin, Inhaler, Simvastatin (as per NOK).	Heparin. Flushed and stored in HTK.
*HMC265	Male	54	C	Head Trauma-MVA	Cigarettes; 1 PPD x 40 yrs.		Fentanyl, insulin, Norcuron, Neosynephrine, Heparin, Albuterol, Dulcolax, Peridex, Docusate, Atrovent, Levaquin, Proteinex, Zantac, Sodium Chloride, Keppra, Tylenol, Midazolam, Ophthalmic Lubricant and Morphine. Flushed and stored in Glut.

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

© DLS

**Donor Information**

Specimen	Gender	Age (years)	Race	Cause of Death	Social History	Medical History	Medication given during Hospitalization
HFC271	Female	49	C	Explosion- Burns	Non-smoker; Non-drinker		Heparin, Regitine, and 25% Albumin.
HMC272	Male	55	C	Head Trauma	1 PPD x 40 yrs. Beer; 6 pck q day x 2 yrs.		Hespan, Unasyn, NSS< Custodial, Lactated Ringers and NaCl. Flushed and stored in HTK.
HMC275	Male	54	C	CVA	Cigarettes- 2 PPD X 20 yrs. Rum and Coke- nightly and Wine coolers > 25 yrs ago; no regular drinking within 15 yrs.	Asthma- treated with inhaler; diagnosed 5 yrs ago. Hypertension; Diagnosed 5 yrs ago; Compliant w/ meds. Diabetes; Diagnosed 5 yrs ago; non-insulin dependent. Pt was treated for Hepatitis in his late teens-type in unk. Blood was found in pt's urine in Feb. 2008. HTN, Cholesterol and Diabetes meds (non insulin), asthma inhaler and nasal spray (as per NOK)	Morphine, Levophed, Neosyneprine, Dopamine, Vasopressin, Kphos, Ancef, solumedrol, and Cleocin, Flushed with UW solution.
*HFC286	Female	52	C	ICH-Stroke	Cigarettes-at least 1 PPK X 35+ yrs. White Russian- 1 every 6 mo. X 35 yrs. Marijuana- Daily X 35+ yrs. Meth-Smoked 1xweek X 10 yrs.	IBS-Diagnosed 30+ yrs ago. Untreated. Ibuprofen and Aspirin (as per NOK)	Insulin, T4, Dopamine, Neosyneprine, Lasix, DDAVP, Cipro, Electrolytes, D5W, K-Phos, and NS (administered in hospital) Flush and storage in UW solution.
*HFC289	Female	65	C	CHI	Cigarettes- 1-2 q d; duration unknown but pt quit since 1989. Wine- 1 glass (since adulthood on a rare basis.)	PT had cervical CA 20 yrs ago. Ieffexor (as reported by NOK.)	Insulin, Phospheryntoin, DDAVP, Hydrocortisone, MGS04, Mannitol, Sodium bicarb, Epinephrine, Fentanyl, Phenylephrine, Levothyroxine, Atomadate, SCCS (DA medication), Versed, Vecuronium, T4, Pitressin, NS, Zosyn and Vasopressin. Insitu: LR/UW and stored in UW solution.
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, Cyrstaloids, Levophed, and Heparin. Epinephrine and atropine were administered. Flushed and stored in UW solution.
HMC297	Male	48	C	Head Trauma-SIGSW	Non-smoker; Non-drinker	Lipitor, Ativan and Seroquel (as per NOK).	Dopamine, Vasopressin, Cefepime, Solumedrol, T4, Narcan, Crystalloids, Albumin, Heparin, Regatine, and Mannitol. Flushed and stored in UW soln.
HFC300	Female	63	C	Anoxia	Beer- 3/day x 30 yrs.	Diabetes; Diagnosed 20yrs ago. Tx w/ Insulin Injections x 15 yrs.	Levophed, NS, Dopamine, Neosyneprine, Heparin, Insulin, Aztreonam, Decadron, Ancef, Nystatin, Erythromycin, Neomycin, Clindamycin, and Magnesium.

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

© DLS

**Donor Information**

Specimen	Gender	Age (years)	Race	Cause of Death	Social History	Medical History	Medication given during Hospitalization
HFC301	Female	45	C	Anoxia	Cigarettes- 1 PPD X >10 yrs. Liquor- Daily. Cocaine- Snorted X 2 yrs. Marijuana- Smoke X 2 yrs. Possibly Crack- X 2 yrs.	Enlarged heart- Diagnosed 2 yrs prior to death. Chest pains- Diagnosed 10 yrs prior to death. Diagnosed with Hypertension 2 years prior to death. Nexium (as per NOK). History of Alcoholism, Benzodiazepine, Methadone, and Opiate abuse.	Piperacillin, Vancomycin, Potassium Chloride, NS Bolus, Albuterol, Zosyn, 0.9 NS, Levophed, PRBCs, NS, Heparin, Mannitol, and Morphine. Flushed and stored in HTK. Epinephrine and Atropine administered at time of CPR
HMC302	Male	49	C	CVA/Stroke-Blunt Injury	Cigarettes- 1 PPD X 30 yrs. Beer- 2-3 / Day X 30 yrs. Marijuana- Smoked as a teen.	Hypertension; Diagnosed 3 yrs ago. Atenolol (as per NOK).	Dopamine, Levophed, Epinephrine, Neosynephrine, Vasopressin, Synthroid, Insulin, Nipride, Hydrocortisone, Zosyn, Lasix, Albumin, RBC, FFP/Plasma, Solumedrol, Mannitol, and Heparin. Flushed and stored in HTK.
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
*HMC310	Male	41	C	Head Trauma-Blunt Injury	Cigarettes- 1/2 PPD since age 15. Beer- 1-2/Weekdays, 9-10/Weekends X 10 years.	COPD- Diagnosed within last 2 yrs. Pneumonia and Bronchitis within last yr. Hypertension; Diagnosed 2 yrs ago. Atenolol, Zyrtec, and Albuterol Inhaler (as per NOK).	Nicardipine, Insulin, Crystalloids, Albuterol, Clindamycin, Gentamycin, Keppra, Rocephin, and Zosyn. Flushed and Stored in HTK.
*HFC313	Female	63	C	Head Trauma-MVA	Wine- 1 Glass/Month.	Possible Basal Cell Carcinoma on Nose. Echinacea and Astragalus (as per NOK)	Dopamine, Vasopressin, Zosyn, KCL, CaCl, Lasix, Insulin, T4, Heparin, Mannitol, and Crystalloids. Flushed and stored in UW solution
*HFC314	Female	71	C	ICH-Stroke	Vodka- 1/2 Pint- Unknown Frequency X 30 yrs.	Pre-Cancerous Mole removed 10 yrs prior to death. Oxycontin, Oxycodone, and Premarin (as per NOK)	PRBC's, FFP, Hespan, Levophed, Vasopressin, Electrolytes, T4, Solumedrol, Sodium Bicarbonate and Labetolol. Flushed and stored in HTK
*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

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

© DLS

**Donor Information**

Specimen	Gender	Age (years)	Race	Cause of Death	Social History	Medical History	Medication given during Hospitalization
*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
*HMC318	Male	53	C	Head Trauma-Blunt Injury	Marijuana; Smoked unknown amount for unknown length of time.	Normal healthy	0.9% NS, Mannitol, Versed, Fentanyl, Levophed, Vasopressin, Pitressin, PRBC's, Magnesium Sulfate, Ancef, Solumedrol, Bicarb, FFP, and Heparin. Flushed and stored in UW solution.
*HFC320	Female	64	C	ICH-Stroke	Cigarettes: < 1 PPD. Quit in 80's. Wine: Quit over 10 yrs prior.	A-Fib: Diagnosed 2 months prior to death. Hypertension: Diagnosed < 5 years prior. Coumadin, Rythmol, Digoxin, Lipitor, Omeprazole and Tramadol (as per NOK)	Levophed, Zemuron, Atropine, Atrovent, Decadron, Desmopressin, Dilantin, K Phos, Nexium, Potassium Phosphate, Crystalloids, Accucheck, D5 1/4 NS, Meropenem, Insulin, KCL, Solumedrol, Magnesium Sulfate, Vitamin K, Lasix, Neosynephrine, Pancuronium, Fentanyl, Heparin and Mannitol. Flushed and stored in Custodial
*HFC321	Female	56	C	ICH-Stroke	Non-smoker; Non-drinker	Arterial Sclerosis: Diagnosed 12 yrs prior. Blood Thinner, Lasix, Pain Meds and Heart Meds (as per NOK)	Epinephrine, Lasix, Neosynephrine, Levophed, Solumedrol, Vasopressin, Ancef, KCL, Erythromycin, Neomycin, Nystatin, Lidocaine, FFP, 0.45 NS, Heparin and Pancuronium. Treated with Vitamin K and FFP.
*HFC323	Female	70	C	CVA-Stroke	Alcohol use.	Asthma: Controlled with Advair and Spiriva. CHF: Diagnosed in Sept 08. Edema: Diagnosed 20+ yrs prior. Mild Hypertension controlled with diet and exercise. Dementia: Diagnosed within the last 2 yrs. Brain Tumor: Diagnosed at age 13. Tumor was removed and patient received radiation X 6 wks. Tumor discovered to be benign. No reoccurrence. Diabetes: Diagnosed 5 yrs prior. Controlled with diet and exercise. GERD: Diagnosed 5 yrs ago. Blood stools Kidney Stones: 20+ yrs prior. Darvacet, Coumadin, Spiriva, Advair, and Nitroglycerine (as per NOK)	evophed, Protonix, Cipro, T4, Solumedrol, Insulin, Lasix, Labetolol, Mannitol, Vasopressin, Digoxin, FFP, PRBC's, 5% Albumin, 1/2 NS, 20 KCL, Heparin, Hextend and CaCl. Flushed with UW, Insulin, Dexamethasone, and Penicillin

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

© DLS

**Donor Information**

Specimen	Gender	Age (years)	Race	Cause of Death	Social History	Medical History	Medication given during Hospitalization
*HMC328	Male	55	C	ICH-Stroke	Whiskey; 4-5 glasses twice a week X 5 yrs.	Vitamins and Aspirin (NOK).	Crystalloids, DDAVP, Lipitor, Norvasc, Vancomycin, 1/2 NS, Dopamine and KCL. Flushed and stored in HTK.
*HFC329	Female	30	C	Anoxia-2 <sup>nd</sup> 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, Acyclovir, Azithromycin, Fexofenadine, Heparin, Hydrochlorothiazide, Dopamine, Lasix, Magnesium Sulfate, metolazone, Metronidazole, Potassium Chloride, Prednisone, Sertraline, Trazadone, Vancomycin and Zosyn. Flushed and stored in HTK.
HMC331	Male	51	C	Anoxia-Cardiovascular	Cigarettes: 1.5 PPD X 30 yrs. Quit in March 2004. Hard Liquor: 2 cocktails/wk X 17 yrs. Marijuana: Smoked 2/ wk X 30 30 yrs.	Shy Drager's Disease; Diagnose August 2008 resulting in Hypotension in addition to some cervical dystonia. Baclofen, Fludrocortisone, Midodrine, Pepcid, Tylenol with Codeine and Vicodin (as per NOK).	Hydrocortisone, Protonix, Depacon, Unasyn, Lovenox, Clonidine, KCL, Lorazepam, K Phos, Dopamine, Lasix, Heparin and Normal Saline. Flushed and stored in UW with Heparin
HMC335	Male	43	C	Head Trauma-MVA	Cigarettes; 1 1/2 PPD since age 15. Beer; six pack and 1/2 since 21. Unsure how long drinking that amount.	Multivitamins (as per NOK).	Neosynephrine, Levophed, T4, Vasopressin, Magsulfate, Calcium Gluconate, Levaquin, Flagyl, Lasix, Insulin, KCL, Kphos, Colloids, Crystalloids, PRBCs, Albumin, NS and Heparin. Flushed and stored in UW soln.
*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.
*HFC338	Female	35	C	ICH-Stroke	Non-smoker; Non-drinker	Certrilene, Ambien and Neurontin (as per NOK)	insulin, T4, Dopamine, Vasopressin, Levothyroxine, Calcium Gluconate, Albuterol, Lasix, Ancef, D5W, PRBCs, Heparin and Mannitol. Flushed and stored in UW.
*HFC339	Female	42	C	Anoxia-CVA	Wine: 10 mL/week.	Cystic Fibrosis, fibroid tumors, and small bowel. CF related Diarrhea. Adek, Multi Vitamins, D plus Calcium, Evening Primrose, Rutin and Ultase (as per NOK)	eftriaxone, Vancomycin, PRBCs, Hespan, NS, Levophed, Kphos, Vasopressin, Insulin, Calcium Gluconate, Dobutamine and Heparin. Flushed and stored in HTK

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

© DLS



**Donor Information**

Specimen	Gender	Age (years)	Race	Cause of Death	Social History	Medical History	Medication given during Hospitalization
*HFC344	Female	50	C	ICH-Stroke	Cigarettes- 1/2 PPD X 35 yrs. Beer- Socially ages 18-35	Allergy related Asthma X 20 yrs. Hypertension- diagnosed 5 yrs prior. Non compliant with meds. Inhaler and HTN meds (as per NOK)	Levophed, Dopamine, Levothyroxine, 1/2 NS, KPhos, Insulin, Solumedrol, D50, 3% Saline, 0.9 % NACL, PRBCs, . 45 NS, D5W, Heparin, T4, Vecuronium and Vasopressors. Flushed and stored in UW solution.
*HFC345	Female	69	C	Anoxia-2 <sup>nd</sup> 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.
*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.
*HFC348	Female	57	C	Cerebrovascular/Stroke	Cigarettes- 1 PPD X 40 yrs.	COPD: Diagnosed 20 yrs prior. Hypertension- Diagnosed 30 yrs prior. Went off meds within last month. Diabetes- diagnosed in her 20s. Tx with meds. Kidney dialysis 3/week X 4 yrs. Pepcid, Dialyvite, Albuterol, and Fosrenal (as per NOK)	Levophed, Crystalloids, Albuterol, Insulin, Kefzol, NS, D5, Dopamine and Heparin. Flushed and Stored in HTK.
HMC349	Male	75	C	Hemorrhagic Bleed	Cigarettes; Quit in early 1990	Angina, CHF. Hypertension; Unknown length of time. Tx with Digoxin. Kidney Stones; early 1960's.	Vitamin K, Fee H2O, Cardizem, Zosyn, Mannitol, Nexium, 1/2 NS, Levophed, Vasopressin, Nipride, Heparin, and Vasopressors. Flushed and stored in UW solution.
HFC350	Female	54	C	ICH-Stroke	Not available.	Multi Vitamins and Calcium (as per NOK). No significant PMH.	Levophed, Levothyroxine, 1/4 NS, Neosynephrine, Lasix, Magnesium Sulfate, Calcium Chloride, Calcium Gluconate, NaHCO3, KPhos, KCL, Vitamin K, Ancef, Solumedrol, Insulin, Heparin, Vasopressors and D50. Flushed and stored in UW solution
*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.

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

© DLS

**Donor Information**

Specimen	Gender	Age (years)	Race	Cause of Death	Social History	Medical History	Medication given during Hospitalization
*HFC352	Female	33	C	ICH-Stroke	Non-smoker; Non-drinker	Seizures since age 2. Became uncontrolled at age 22. Pt is quadriplegic and confined to wheelchair with impulse control disorder. Colace, Cerovite Liquid, Paroxetine, Hydrochloride, Dilantin, Keppra, Topamax, Levothyroxine Sodium, Klor, Milk of Magnesia and Multi Vitamins (as per NOK)	Levophed, Insulin, Crystalloids, Ancef, Calcium Gluconate, DDAVP, Magnesium Sulfate, Potassium Acetate, Potassium Chloride, Vancomycin and Zosyn. Flushed and stored in HTK.
*HFC356	Female	42	C	ICH-Stroke	Cigarettes- 1PPD X 3 yrs. Gin and Tonic- 1/day X 18 yrs.	Hypertension- Diagnosed a few yrs. Prior. Neurological Disease- At age 22 pt had surgery and was given epilepsy med. Condition resolved with no issues since. Tetracycline, Daily Vitamins, and Fish Oil (as per NOK)	Phenylephrine, Vasopressin, Levophed, Dopamine, Neosynephrine, T4, Albuterol, Calcium Chloride, Dextrose, Insulin, Lasix, Sodium Bicarbonate, NS, KCL, Solumedrol, Cipro, Dobutamine, Heparin and Mannitol. Flushed and stored in HTK.
HFC358	Female	52	C	ICH-Stroke	Non-smoker; Non-drinker	Hypertension- Diagnosed 31 yrs prior. Atenolol, Allopurinol, and albuterol (as per NOK)	Levophed, Vasopressin, Labetolol, KCL, Solumedrol, Zosyn, Combivent, Insulin, Lasix, T4, Lopressor, Magnesium Sulfate, KPhos, Crystalloids, Heparin and Solumedrol. 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.
*HMC360	Male	55	C	ICH-Stroke	Cigarettes: 2 1/2 PPD X 35 yrs. Vodka and Beer: 1-2 drinks/ month X 20 yrs.	Hypertension; Diagnosed 1 yr ago. Diabetes; Diagnosed 4 months ago. GERD; Diagnosed 8 yrs ago. Prilosec, Metformin, ASA, Plavix, Zocor, Lisinopril, and Glimepiride (as per NOK).	Cardene, Levophed, 1/2 NS, Ancef and Heparin. Flushed and stored in HTK.
*HFC363	Female	65	C	Cerebrovascular/Stroke	Non-smoker; Non-drinker	Open heart surgery 7 yrs prior. Insulin dependent Diabetes X 30 yrs. GERD. ASA, Metformin, Plavix, Metoprolol, Reglan, Trazadone, Synthroid, Lasix, Prilosec, Welbutrin, Cozaar, Cymbalta, K-Dur, Isosorbide, Evista and Motrin (as per NOK)	Levophed, Neosynephrine, Dopamine, Solumedrol, T4, Insulin, Arginine, Zosyn and Vancomycin (administered in hospital) Flushed and stored in HTK.
*HFC365	Female	42	C	Anoxia- 2 <sup>nd</sup> 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.

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

© DLS

**Donor Information**

Specimen	Gender	Age (years)	Race	Cause of Death	Social History	Medical History	Medication given during Hospitalization
*HFC368	Female	64	C	Head Trauma-MVA	Margarita- Few times a year.	Hypertension-Diagnosed 20 yrs. Prior. BP Meds and Calcium (as per NOK)	Neosynephrine, Vasopressin, Dopamine, Levophed, Insulin, Ancef, Zosyn, Nexium, Cerebryx, Crystalloids, Colloids, T4 and Heparin. Flushed and stored in HTK.
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.
*HFC377	Female	54	C	CVA-ICH	Cigarettes- 1 PPD X 30 yrs. Quit in Jan 2005.	Bypass surgery- 1999. Hypertension- Diagnosed 8 yrs prior. Nitro (as per NOK)	Neosynephrine, Zosyn, Solumedrol, Albuterol, T4, Insulin, Dopamine, Adenosine, Bicarb, 5% Albumin, ICA, Levophed, FFP, PRBCs, Crystalloids and Heparin (administered in hospital) Flushing Solution: UW. Storage Solution: UW.
*HFC379	Female	37	C	ICH-Stroke	Cigarettes; 2-3 PPD x 10-15 yrs. Vodka/wine coolers- on occasion. Marijuana-experimented with in HS.	Hypertension; diagnosed in early 20's. Xanax, BP Meds, and Multi Vitamin (as per NOK).	Dopamine, Levophed, Amiodarone, Heparin, Mannitol, Vasopressors and Crystalloids. Flushed and stored in UW solution.
*HFC380	Female	45	C	Anoxia-2 <sup>nd</sup> to Respiratory Failure	Non-smoker; Non-drinker	Severe Asthma >30 yrs. Has been hospitalized over 10 times for Asthma related reasons.	Levophed, Fentanyl, Cardene, Ativan, Crystalloids, Vasopressin, Zemuron, Combivent, Fragmin, Insulin, Lasix, Levothyroxine, Mannitol, Narcan, Solumedrol, Zosyn, Zyvox, NS, Pitressin, T4, Heparin and Vasopressin. Flushed and stored in HTK.
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.
*HFC385	Female	63	C	ICH-Stroke	Cigarettes- unknown amount or duration. Quit 40 yrs ago.	Coronary artery bypass graph- 6 vessels in 1995. Hypertension- Diagnosed 20 yrs. ago. Diabetes- diagnosed 7 yrs. Prior. Dialysis- 7 yrs ago.	Levothyroxine, Solumedrol, Insulin, Neosynephrine, Ancef, Dopamine, Keppra, Dextrose, Crystalloids, Heparin, T4 and Vecuronium. Flushed and stored in UW solution.

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

© DLS

**Donor Information**

Specimen	Gender	Age (years)	Race	Cause of Death	Social History	Medical History	Medication given during Hospitalization
HFC388	Female	67	C	Head Trauma- 2 <sup>nd</sup> to fall	Wine- Rarely once a month	Breast Cancer- Pt had mastectomy and received Chemo and radiation 15 yrs. Prior. Cancer free 5 yrs.	Levophed, Dopamine, Vasopressin, Levothyroxine, Dobutamine, Magnesium Sulfate, Adenosine, Hydrocortisone, Heparin, Mannitol, Lasix, Solumedrol and Crystalloids. Flushed and stored in HTK.
*HFC430	Female	50	C	Anoxia-2 <sup>nd</sup> to Cardiovascular	Cigarettes- 2 PPD X 25 yrs. Vodka & Mixed Drinks- Daily X 25 yrs. Cocaine & Marijuana- 6 yrs ago.	Asthma- Developed as an adult. Hypertension- Diagnosed 10 yrs prior. Elevated liver enzymes upon admit. HCTZ (as per NOK)	Propofol, Unasyn, D5 1/2 NS, KCL, Metoprolol, Heparin and Thorazine. Flushed and stored in HTK.
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.
*HFC448	Female	71	C	ICH-Stroke	Non-smoker; Non-drinker	Hypertension- Diagnosed 2 months prior. Diabetes- Diagnosed 2 months prior: Insulin dependent. Kidney stones- 30 yrs prior.	Dopamine, Neosynphrine, I.V. Fluids, Insulin and PRBCs. Flushed and stored in UW solution.
*HFC449	Female	57	C	ICH-Stroke	Non-smoker; Non-drinker	Hypertension- Diagnosed 3 yrs prior. Compliant with meds. Recent Seizure upon admit. Type II Diabetes- Diagnosed 6 yrs prior. Compliant with meds. Zydis, Lovastatin, Acetaminophen, Aspirin, Bupropion, Dicyclomine, Escitalopram, Flurazepam, Gabapentin, Lisinopril, Metformin, Ziprasidone (as per NOK).	NS, Dopamine, Levaquin, Clindamycin, Insulin, Levothyroxine, Narcan, Norcuron, Pitressin, Potassium Chloride, Sodium Bicarbonate, Solumedrol, Water, Albuterol, Atrovent, Magnesium.
HFC450	Female	45	C	CVA	1 bottle per wk for the last 10 yrs.	Pt's medical hx significant for migraine.	Dobutamine, Neosynphrine, Levophed, Milrinone, Heparin and Amicar. Flushed and stored in SPS-1
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.
*HFC456	Female	74	C	ICH-Stroke	Cigarettes- Quit 45 yrs prior. Wine, Liquor, & mixed drinks- Very Seldom, Since becoming an adult.	Aortic valve replacement 2000. CABG. Pacemaker implant 1999. Hypertension- Diagnosed 20 yrs prior. Dementia/Alzheimer's- Diagnosed 5 yrs prior. Type I Diabetes- Diagnosed 30 yrs ago- Tx with Insulin Injections. Aricept, Lantus, Lovastatin, Namemdia, Metoprolol. Ramipril, and Warfarin (as per NOK).	FFP, Dopamine, Heparin, and Regitine. Flushed and stored in UW solution.

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

© DLS

**Donor Information**

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

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

© DLS

**Donor Information**

Specimen	Gender	Age (years)	Race	Cause of Death	Social History	Medical History	Medication given during Hospitalization
HFC482	Female	63	C	Cardiovascular 2 <sup>nd</sup> ICH	Cigarettes: 2 PPD since 15 yr old. Tequila: 1 time per week X 5 yrs, amount unknown. Pt quit drinking in 2002	Uterine cancer at age 20 and breast biopsy 3 yrs prior which was benign. Bloody stools from 1997-2002, only when drinking. Pt received several colonoscopies with no malignancy found. Lortab, pain medication, estrogen, prescribed but non-compliant (as per NOK). Pt was allergic to Barium and penicillin (as per NOK)	Heparin, Protamine, NaCHO <sub>3</sub> , Epinephrine, KCL, K Phos., Levophed, Neosynephrine, 1/2 NS, 1/4 NS, LR, Haspan, Pitressin, Methylprednisolone, K+, FFP, PRBCs, Platelets, Tylenol, Magnesium, Solumedrol, Insulin, Levothyroxine, Avelox, Sodium Bicarbonate, Dobutamine, Narcan, Vasopressin, Crystalloids, VitaminK, Calcium Chloride, Albumin, Norepinephrine, Mannitol, Lasix, Diuril, Magnesium Sulfate, Ephedrinem and Phenylephrine. Flushed with HTK and stored in UW
*HFC483	Female	76	C	ICH-Stroke	Cigarettes: at least 1 PPD X 60+ yrs. Beer: at least 3 per day X 60 yrs.	Diagnosed with Hypertension 8-9 yrs prior. Med compliant. Unknown blood thinner ( as per NOK)	Phenylephrine, Cefazolin, NS, Labetolol, Glucose, Heparin, Ephedrine, Ondanstron, Nicardipine, Folic Acid, multi vitamin, Pantaprazole, Etomidate, Succinylcholine, Arovastatin, Docusate, Hextend, Insulin, T4, Neosynephrine and Vasopressin. Flushed and stored in HTK.
*HFC485	Female	56	C	ICH-Stroke	Filtered Cigarettes- 1 Pack per week X 5 yrs. Wine - 1 glass & only during holidays X 30 yrs.	Hypertensive. 4 Blood pressure medications - exact names unknown (as reported by NOK)	Levophed, Neosynephrine, Gentamycin, Mannitol, Daptomycin, Sodium Bicarbonate, KPfos, and Magnesium. Flushed and stored in HTK.
*HFC488	Female	43	C	Anoxia-2 <sup>nd</sup> to Cardiac Arrest	Cigarettes: 4-5 ciggs X 3 yrs. Beer: 1-2 rarely X 25 yrs.	Normal healthy	Fentanyl, Sodium BiPhos, Glutamine, Protonix, Lovenox, Oxycodone, Ampicillin, Chloraseptic, Vancomycin, Zosyn, Tylenol, Motrin, Sodium Chloride, Insulin, RBCs, FFP, Dopamine and Levophed. Flushed and stored in LR/UW
HMC489	Male	82	C	CVA-Stroke	Non-smoker; Non-drinker	Hypertension: Diagnosed 10-15 yrs ago. Type II Diabetes: Diagnosed last October. Bleeding Ulcers: 1953- No reoccurrence. Avodor, Glemepeside, Simvastin, Enapril/HCTZ and Lubiprostone (as per NOK).	T4, Neosynephrine, Crystalloids, Ancef, Heparin, Mannitol, and Lasix. Flushed and stored in UW solution.

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

© DLS

**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.
*HFC499	Female	28	C	Head Trauma-2 <sup>nd</sup> to MVA	Cigarettes: 2.5 PPD X 13 yrs. Beer: Bottle a day on weekends. Marijuana: Smoked every 3 months X 10 yrs. Last used a week prior.	Mild Bronchitis- 2-3 times a year. Treated with Duoneb. Pre Clampsia- With both pregnancies. Last time 6 months. Fish Oil (as reported by NOK)	V Dextrose, Insulin, PRBCs,
*HMC500	Male	54	C	CVA-ICH	Beer: Occasionally 1-2 times a week since 18 yrs old.	CVA- 3 months ago requiring 1 wk hospital stay. CVA: 7 yrs ago. Lexapro, Lipitor, ASA, Altace, Coumadin, and Alovert (as per NOK).	Crystalloids, Colloids, Cardene, Vancomycin, Platelets, and Heparin. Flushed and stored in UW solution.
*HFC506	Female	72	C	Anoxia-2 <sup>nd</sup> to Cardiovascular	Social alcohol at holidays X adult life.	History of Asthma- Treated with Advair. Hypertension- Diagnosed > 40 yrs ago. Treated with Norvasc. History of Gait changes- Vertigo. Type II Diabetes X 15 yrs. Insulin dependent X 15 yrs. Irritable bowel syndrome. Kidney Failure X 4 yrs. Kidney Stones over last couple years. Treated with Kidney Dialysis in 2001. Frequent bladder infections & UTIs. Norvasc, Advair, Acetaminophen, Aspirin, Florajen, Gemfibrozil, Humalog, Lantus, Percocet, Prednisone, Prograf, Bactrim and Toprol as per NOK).	IV Fluids, Methylprednisone, Hydrocortisone, Insulin, Vasopressin, Dopamine, Levophed, Bicarb, Heparin, Calcium, Invanz, Lasix, Albumin, Magnesium, Kphos, KCL, Azithromycin, Vancomycin, Meropenum, Pitressin, CaCl, LR, 0.9 NS, and Hespan. Flushed and stored in UW solution.

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

© DLS



**Donor Information**

Specimen	Gender	Age (years)	Race	Cause of Death	Social History	Medical History	Medication given during Hospitalization
*HFC507	Female	61	C	ICH-Stroke	Cigarettes- Smoked < 1 Pack per yr socially. Alcohol use very infrequently: once a year.	History of CHF, Pacemaker, and A-Fib. Hypertension- Diagnosed 35 years- Compliant with meds. Diabetes- Type II X 2-3 yrs. Took oral meds & controlled with diet. Gabapentin, Lisinopril, Januvia, Metoprolol, Metolazone, Hydralazine, Bumetanide, Klor Con, Isosorbide, Coumadin, ASA, and Vege Lax (as per NOK).	Dopamine, IV Fluids, Lasix, Bumex, T4, FFP, Bicarb, Neosynephrine, Levophed, Dobutamine, Epinephrine, and Heparin. Flushed and stored in UW solution.
*HMC510	Male	22	C	Head Trauma-2 <sup>nd</sup> 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.
*HMC513	Male	68	C	ICH-Stroke	Cigarettes: 1 PPD since age 17. Quit in 1986. Beer: 1 a week socially since a teenager.	COPD: Treated with meds due to heart problems. CAD/CHF since 1987. CABG X 2. Stents. History of LAVD. Hypertension: Diagnosed in 1987. Diabetes in 1987. Treated with Novolog and Lantis. Insulin dependent X 20 yrs. Meds for HTN, Diabetes, Depression, Heart, Multivitamin, Magnesium, Folic Acid and Protonics (as per NOK). History of LVAD and Von Willebrand Disease.	Propofol, Heparin, Coumadin, Insulin, Levophed, amitriptyline, Aspirin, Carvediol, Cetrizine, Divalproex, Docusate, Folic Acid, Hydrazaline, Hydrocodone, Insulin Glargine, Insulin Lispro, Isosorbide, Lisinopril, Magnesium Oxide, Magnesium Sulphate, 25% Mannitol, Methimazole, Morphine, Multivitamin, Naloxone, Pantoprazole, Potassium Chloride, Prochlorperazine, Rosuvastatin, Topiramate, Venlafaxine, Warfarin Sodium, Spironolactone, FFP, 5% Albumin, NS, Pitressin, Narcan, Solumedrol, KCL, Kphos, Ancef, Heparin, and Hespan. Flushed and stored in UW solution.
*HFC518	Female	42	C	CVA 2 <sup>nd</sup> ICH	Non-smoker; Non-drinker	Donor was prescribed Paxil; reason not mentioned in NOK questionnaire. Donor took antidepressant Paxil, Ambien as needed and not sure if still taking cholesterol medication (as per NOK).	Neosynephrine, Levophed, Insulin, IVPB, Vasopressin, 1/2 NS, D5 1/4 NS/20kcl, 1/2 NS/bicarb, KCL, Lasix, DDAVP, Albuterol, Unasyn, Levothyroxine, Cefalozin, Solumedrol, Levaquin, Narcan, Heparin, Crystalloids, Epinephrine, Norepinephrine, and Dopamine, Flushed and stored in SPS.

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

© DLS



**Donor Information**

Specimen	Gender	Age (years)	Race	Cause of Death	Social History	Medical History	Medication given during Hospitalization
*HFC519	Female	54	C	CVA 2 <sup>nd</sup> ICH	Non-smoker; Non-drinker	Donor had no known cancer, however, suffered with benign ovarian cysts which led to a hysterectomy. Donors chart shows "insulin resistance" and DM type 2. NOK states none in medical/social history. Donors medical/social history states NO liver disease, however, interviewer mentions in notes further complications of liver biopsy in 2009. Donor was diagnosed with Celiac Disease in 2009. Lisinopril, Zanaflex, Remeron, Urodiol, Ritalin, Klor-con, Klonopin, Levothyroxine, Lasix, Trazadone, and Nitrofurantoin (as prescribed by physician).	Levophed, Bicarb, Xigris, Solumedrol, insulin, Vecuronium, Zosyn, Zithromax Succinylchloride, Etomidate, Narcan, Rocephin, Heparin and Crystalloids. Flushed and stored in UW solution.
HMC521	Male	25	C	SIGSW	Abuse rehab 7 years ago; alcohol, cocaine (inhaled). Positive for alcohol on admission.	Herpes.	Dopamine, D5, KCl, potassium, Ancef, DDAVP, Hespan.
HFC529	Female	64	C	CVA 2 <sup>nd</sup> ICH	Pt smoked for 3-4 yrs during college but quit after graduating and has not smoked since. Alcohol; Seldom and on special occasions.	Pituitary tumor; Diagnosed 15-20 yrs ago. It was found to be benign and checked annually. Pt took Vitamins and a calcium supplement daily (as per NOK).	Pitressin, D5, Nicardipine, Insulin, Narcan, Hydralazine, NS, Ancef, Norcuron, Atrovent, Albuterol, Humulin, Cardizem, Vecuronium, Neosynephrine, Solumedrol, Insulin, Vasopressin, Heparin, Antihypertensives. Flushed and stored in HTK.
*HMC530	Male	51	C	Anoxia-2 <sup>nd</sup> to Cardiovascular	Mixed Drink; 1 every 6 months.	A-Fib; Diagnosed 1 yr prior. Leaky Valve Diagnosed 1 yr prior. Hypertension; Diagnosed 1 yr prior. Diabetes; undiagnosed (NOK claims diet controlled "sugar problem.") Cardizem, Digoxin, Aspirin, Lasix, Klorcon, Metoprolol, and Pantopzoic (as per NOK).	Levophed, Crystalloids, Colloids, Amiodarone, Vancomycin, Ancef, Platelets, and Heparin. Flushed and stored in UW solution.
*HMC554	Male	46	C	Anoxic injury	Normal.	IDDM since age 12. Appendectomy 25 years ago. MI 1992. Ulcers. Bilat AVF. Kidney dialysis since 3/92, received kidney transplant 8/13/95. Vaccinated for Hepatitis B 5/92. Atonally, Basal gel, insulin, Pepcid, Quinine, Coumadin, Niferex, Flourinef, ISMO.	Dobutamine, insulin, NaCl, Lidocaine, NT6, Nepinephrine, Heparin, Dopamine, crystalloids, colloids, blood products, Acyclovir, Azathioprine, Diphenhydramine HCl, Lasix, Solumedrol, Sulfamethoxazole, Bactrim, Metoclopramide, Antithymocytic globulin, Cleforan.
*HFC580	Female	64	C	Cerebrovascular intracranial hemorrhage	Past heavy smoking and drug abuse, requiring hospitalization '95. Placed in nursing home where did well with recovery.	Had appendix surgery as a child, broken ribs 15-18 yrs ago. Hospitalized in past two years for Jaundice.	Dopamine, Nipride, Blood Products, Crystalloids

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

© DLS

**Donor Information**

Specimen	Gender	Age (years)	Race	Cause of Death	Social History	Medical History	Medication given during Hospitalization
*HMC592	Male	56	C	Anoxia, carbon monoxide poisoning	Social alcohol. Military service x20 years, in Vietnam twice (not currently active military).	Melanoma, removed from face in last year. Acute arthritis. STDs in past. Achilles tendon repair 1987, 1989. Erythromycin, Robitussin for upper respiratory infection. Possible antidepressant.	Heparin, Solumedrol, Lidocaine, Dopamine, Dobutamine, Hespan, NaBicarb, crystalloids.
*HMC594	Male	70	C	CVA	Social alcohol use.	Craniotomy as child. Arthritis in neck, back. Disc disease. Non-spontaneous arrhythmia. Frequent GI ulcers. Zantac, Inderal.	Dopamine, Nipride, Neosynephrine.
HMC596	Male	44	C	SAH s/p aneurysm	Social alcohol. Marijuana.	Mumps as child. Shortness of breath with activity. Borderline mental retardation. Klinefelter's disease. Hormone replacement x2-3 months 10 years ago. Hypertension, dx 2-3 weeks prior to admission. Medizone for nausea/vertigo 3 days prior to admission.	Lasix, Dilantin, Dobutamine, Decadron, Vasopressin, Pitressin, Nimodysine, T4, Solumedrol, Ancef, Oracillin, Alupent, Tylenol, Ampicillin, Gent, Clindamycin, Lidocaine, Thio, Pavulon, Succs, Ativan, Timintin, Vancomycin, CaGluconate, KCl, D5, crystalloids.
HFC603	Female	65	C	Intracranial hemorrhage	Smoker 3ppd x 40 yrs	Hypertension x4 years. Echo 1995. 2 cardiac cath, last one in 1992. EKG 1995. Stress test 1983, 1987. Appendectomy 1948. Gallbladder 1951. Scoliosis. Hysterectomy 1977. Colitis, diverticulitis. Bowel resection 1979. Breast biopsy 1995, results negative. Degenerative, osteo arthritis. Hepatitis 1996. Skin graft on hand. Elective spinal fusion surgery on admission. Beginning stage emphysema. Zoloft, Vasotec, Losol, Premarin, Ativan, Elavil, Darvocet. Food allergies, allergic to codeine. Serologies indicate previous hepatitis exposure or immunity (note: PCR test is negative).	Neosynephrine, Norepinephrine, Mannitol, Oxacillin, Pitressin.
*HMC605	Male	59	C	Cerebrovascular Accident / hemorrhage	Quit drinking 1992; prior, alcohol abuse 30-40 years.	Asthma, treated with inhalers. COPD. Angina, CABG 1987. Hypertension x1987. Atenol, Nifedipineim, ASA, Isordil (recently stopped).	Heparin, Dopamine, Esmolol, Nipride, TPA, nitroglycerine, Diltizem.
HMC608	Male	45	C	Subarachnoid hemorrhage s/p fall from horse	Heavy drinker, more so last few years; 12 beers/day, +liquor. Marijuana many years ago. Possible jail time over 19 years ago for speeding.	Skin graft 24 years ago. Broke foot 8-10 years ago. Occasional heartburn; Roloids, Tagamet. Vitamins, Geritol. Craniotomy on admission.	Heparin, Mannitol, Vasopressin, Dopamine, Pavulon, Levophed, Tylenol, MSO4, albumin, crystalloids.
*HMC627	Male	44	C	Anoxia, cardiorespiratory arrest	Normal.	Acid peptic disease, possibly pancreatitis. Questionable GI bleeding 10-15 years ago. Mentally handicapped, cerebral palsy, lived in group home with Level 4 care. State hospital from ages 16 to 21. Tagamet, Clozaril, Mylanta. Possible allergy to Thorazine.	Decadron, Rocephin, Ancef, Cardizem, Dopamine, Epinephrine, Atropine, Lidocaine, NaHCO3, crystalloids, sodium bicarb.

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

© DLS

**Donor Information**

Specimen	Gender	Age (years)	Race	Cause of Death	Social History	Medical History	Medication given during Hospitalization
HMC633	Male	26	C	Self inflicted gun shot wound	Marijuana, heavy drinker.	Cut fingers/hand in last two years.	Dopamine, Dobutamine, Pitressin, Levothyroxine, Solumedrol, D50, insulin, Zinicef, Lasix, crystalloids, colloids.
HMC634	Male	57	C	Subarachnoid hemorrhage	Whiskey.	Heart disease >10 years. Arthritis, stiff joints. Hernia operation 13 years ago. Arthroscopic knee surgery 5 years ago. Indoral, HCTZ, aspirin. Amoxicillin x2 weeks.	Heparin, Epinephrine, Zemuron, Dopamine, Hespan, Piperacillin, Cipro, Diflucan, crystalloids, colloids.
*HF644	Female	76	C	Anoxia, CVA	1-2 drinks per week socially.	D and C 1960. History of herpes. Recent hacking cough. Recent slow heartbeat. Arthritis, some back pain. Pseudoephedrine. IIAM notes state non-malignant tumor discovered on uterus during organ recovery; no medical forms to confirm.	Dopamine, Atropine, Epinephrine, crystalloids.
*HMC657	Male	64	C	Intracranial Bleed	Heavy alcohol use in past, cut down in last year.	Hypertension. Thyroid problems. Asbestos exposure 30 years ago. Blood in urine 2 weeks ago; renal carcinoma noted on kidney biopsy. Synthroid, Glynase, Ceflex.	Nipride, Deiprivan, Decadron, Pepcid, Dilantin, Nitro paste, Nimodipine, Ancef, KCl, Mg
*HMC662	Male	18	C	Closed head injury s/p motor vehicle accident	Social alcohol use.	Lumbar laminectomy 4 years ago. Several bouts of pneumonia, diagnosed with pleurisy several weeks prior to admission.	Heparin, Mannitol, Rocephin, Pepcid, Timentin, Lasix, Solumedrol, Motrin, Tylenol, Ancef, Neosynephrine, Dopamine, Neomycin, Erythromycin, Nystatin, albumin, blood products, crystalloids, Unasyn, Fortaz, Norcuron.
HFC664	Female	50	C	Subarachnoid Hemorrhage	Smoker, drinker (10 beers/wk x 15 yrs), marijuana a few times. Tattoo left chest, professionally done.	Heart murmur, Tachycardia 20 yrs ago, complained of palpitations from time to time. Gall bladder surgery 10 yrs ago. Knee joints popped all the time.	Aqueous, Pitressin, Ancef
HMC678	Male	20	C	Head trauma s/p Motor vehicle accident.	Occasional 1 pd, Social alcohol. History of cocaine, marijuana (no IV drugs); clean for months. Tattoos, professional. Jail, 4-5 days, 1 year ago for B&E.	Juvenile diabetes, oral insulin x4 years. Hernia repair as infant. Broke right knee 3 years ago. Orchiectomy. Tonsils, adenoids. Craniotomy and splenectomy on admission.	Heparin, Thyroxine, Kefzol, Dopamine, Levophed, Lasix, Potassium, Dilantin, Pitressin, T3, Neosynephrine, insulin, albumin, crystalloids, blood products.
*HMC686	Male	56	C	Intracranial hemorrhage s/p motor cycle accident	Social alcohol. Hospitalized for depression, OCD 2/97.	Varicose veins 1969. Hemorrhoid 1971. Depression, OCD. Possibly borderline diabetic.	Dopamine, Levophed, Lidocaine, D5, albumin, crystalloids, colloids.

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

© DLS

**Donor Information**

Specimen	Gender	Age (years)	Race	Cause of Death	Social History	Medical History	Medication given during Hospitalization
*HMC693	Male	72	C	Anoxia	Social alcohol use.	Asthma, seasonal hay fever. Hypertension x20 years. Arthritis in back. Aortic aneurysm 6 years ago. Cholecystectomy. Gastrectomy. Lower leg cellulitis 1 year ago, treated with antibiotics. Arthroscopy for cardiac arrest. HCTZ, Tenormin, Vasotec. Thoracic aneurysm repaired, CABG on admission.	Heparin, Mannitol, Versed, Fentanyl, Dopamine, Lidocaine, Dobutamine, Epinephrine, Labetolol, Atenolol, Enalapril, Tylenol 6, Lorazepam, aspirin, Neosynephrine, Pancuronium, Protamine, sodium bicarb, Cefazolin, Bemetanide, CaChloride, Lidocaine, Temazepam, sodium chloride, insulin, crystalloids, colloids.
*HFC699	Female	60	C	Cerebrovascular Accident, Intracranial hemorrhage	Normal.	Flu-like symptoms. Brain aneurysm dx in last 6 weeks. Rheumatic arthritis x16 years. Hysterectomy. Appendectomy. Tonsillectomy. Aspirin, Methotrexate, estrogen patch.	Heparin, Mannitol, Dopamine, Pitressin, D5, Nipride, Nimopridine, NTG, Pavulon, Epinephrine, Decadron, Fosphenytoin, Esmolol, insulin, Isoflurane, Ancef, KCl, CaCl, Pepcid, crystalloids.
HFC707	Female	60	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
*HMC712	Male	45	C	Head Trauma	Smoker, Infrequent vodka drinker	Cluster migraines. Elevated blood pressure 2nd to anxiety. Substance abuse problem last 10 yrs for Tylenol w/codeine, hydrocodine, Percocet off/on. Daily meds Anacin, Aleve, Prozac, Catapress, Tylenol w/codeine, Percocet	Dopamine, Mannitol, Decdron, Zantac, Pitressin, Succinylcholine
HFC716	Female	73	C	Cerebrovascular Accident	Infrequent Smoker, Social drinker	Glaucoma, Spastic colon w/Esophagitis, gastritis. Fibromyoglia. HTN since 1985. Prior surgeries appendix removal, Hysterectomy.	Dopamine, Nipride, Vitamin K, Magnesium, Dobutamine, Zantac, Pitressin, Solumedrol, MgSO4, Hespal, Neosyndrine, Phosphate, Calcium, Heparin
HFC725	Female	70	C	Stroke, ICB	2 glasses of wine before dinner, 1 Scotch daily.	Hypertension x15 years. Cholecystectomy 1971. Ankle fracture 1989. Normadyne, Antivert. Allergic to sulfa drugs.	Heparin, Dopamine, Levaquin, D5, Epinephrine, Atropine, albumin, crystalloids, colloids
*HFC726	Female	71	C	ICH	3 or 4 drinks a week.	History of arteriovenous malformations. Prior ICBs 1965, 1967, 1976. Knee replacement 1997. Mitral valve replacement 1995. Hemorrhoids. Skin cancer on nose 5 or 6 years ago; malignant, not melanoma. Cataract surgery 2-3 years ago. Coumadin, Detrol, Celebrex.	Heparin, Dopamine, KCl, Ancef, insulin, D5, Vit K, blood products.

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

© DLS

**Donor Information**

Specimen	Gender	Age (years)	Race	Cause of Death	Social History	Medical History	Medication given during Hospitalization
*HMC728	Male	78	C	ICH	Social alcohol use.	Hypertension, recently diagnosed. Arthroscopic knee surgery 1982. Polyps removed 1999. Basal tumor on nose 30 years ago. Bypass surgery 1995. Lopressor, Digoxin, Coumadin, Metoprolol, Lipitor, Maxide. Possibly vaccinated for Hep B during WWII in service.	Heparin, Dopamine, Lidocaine, Vecuronium, Ancef, CaGluc, KCl, crystalloids, blood products.
*HFC732	Female	41	C	Intracranial Hemorrhage	Clean 2 years. Quit drinking 2 years ago, prior fifth of liquor, 6 pack beer daily. Tattoo (nonprofessional). Marijuana x8 years, quit 2 years ago. Speed x28 years, quit 2 years ago. Boyfriend jailed overnight, 2 years ago, drug charges.	Borderline diabetic >20 years. Bulimic dx 2 years ago, likely x5 years. Chemical exposure at work (furniture company) 8 years ago. Asthma meds for less than 1 year subsequent to chem exposure. Multiple broken >20 years ago. Uterine cancer 13-15 years ago, partial hysterectomy. Seizures, unknown etiology >20 years, worsening last 5 years, associated memory loss. Dilantin, Tegretal.	Heparin, Mannitol, Dopamine, Dobutamine, Dilantin, Neosynephrine, T3, T4, Rocuronium, Synthroid, Ancef, CaCl, albumin, crystalloids.
*HMC733	Male	30	C	Head injury s/p motor vehicle accident	Beer, 6 pack at a time, not daily, x13 years. Drinking moonshine prior to admission. Cocaine, snorted, marijuana. Jail, 5/00 sentenced to 50 days, 1 1/2 years ago served 15 months, probation violation both times. Tattoos.	Metal plate in skull, from skull fracture 11 months previous. Valium. Some GI bleeding x1 year, doctor said ETOH related. Broke wrist 15 years ago.	Dopamine, Vasopressin, blood products, crystalloids, colloids, albumin, Levothyroxin, KCl, Ancef, Bumex, Lasix, tetanus, MgSO4, CaGI, HCO3, Solucortef, Mannitol, Nimbox
*HFC736	Female	67	C	Intracranial Hemorrhage	Ex-smoker, Non-drinker	Hysterectomy Glaucoma Daily meds drops for glaucoma.	Ancef, Dopamine, DDAVP, Heparin, Mannitol, Lasix, Solucortef
*HFC737	Female	77	C	Cerebrovascular Accident	Non-smoker, occasional glass of wine or beer.	Broken foot 1958. Choliapertomy 30 yrs ago. Daily meds Denvasc, Digoxin, Ecolrin, Premarin	decadron, Pepcid, Lasix
HMC753	Male	55	C	Intracranial hemorrhage s/p cerebral edema	2-3 wine/liquor per day.	Acid reflux, treated with Pepcid. Broken leg, dislocated shoulder, deviated septum. Possible active infection at TOD - meningitis/encephalitis.	Dopamine, Rocephen, Ampicillin, Vancomycin, Acyclovir, Decadron, Ativan, blood products, Heparin, Vecuronium, morphine, Diprovan, Lasix, Propofol, Pepcid, Nipride
HMC760	Male	56	C	Head trauma s/p motor cycle accident	Normal.	Head surgery 1985. Seasonal asthma, used inhaler. Biopsy notes low density lesion on left lobe of liver, possible hemangioma.	Heparin, Dopamine, crystalloids, colloids
*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

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

© DLS

**Donor Information**

Specimen	Gender	Age (years)	Race	Cause of Death	Social History	Medical History	Medication given during Hospitalization
*HMC767	Male	46	C	Cerebrovascular Accident, Intracranial hemorrhage	Quit drinking 20 years ago. Occasional marijuana, quit 15 years ago. Tattoos, self-administered >26 years ago. Pierced ear, self-administered, 5 years ago.	Allergic to dog hair. Torn cartilage in knees from football. Broke hand 5-6 years ago. Hypertension, controlled by meds (unknown). Excedrin; 2 years ago took herbal supplement super blue green algae.	Propofol, Catapres, Labetalol, Mannitol, Dilantin, Unasyn, Solumedrol, KCl, D5W, Nipride, insulin, Heparin, crystalloids.
HMC775	Male	76	C	CVA-ICH	Normal.	Persistent cough, shortness of breath secondary to cardiac insufficiency. CHF, A fib. CABG, valve replacement 1/01. Cardiac cath 12/00. Anticoagulation therapy. Yellow jaundice during WWII. Nonmalignant polyp removed 8-10 years ago. Charcot Foot syndrome. Hypertension, controlled by meds, >8 years. Hypothyroidism. Macular degeneration. Bladder problems. Prostate surgery, hemorrhoidectomy 1991.. OTC anti inflammatories for arthritis. Asbestos exposure in work building. Recent prescription for Amoxicillin for dental work. Ditropan, Lasix, Pacerone, Diovan, Coumadin, Lanoxin, Synthroid.	Heparin, Norcuron, Lidocaine, Dopamine, Clindamycin, DDAVP, KCl, crystalloids.
*HMC777	Male	35	C	Anoxia, ruptured aorta s/p Motor Vehicle Accident	Beer 3x week, >5 bottles each time. Marijuana, cocaine, steroids, all 15 years ago. Worked for septic company - exposure to medical waste. Tattoo.	Depression x1 month. Effexor. Possibly vaccinated for Hep B.	Heparin, Esmolol, Labetalol, Nitroglycerine, Ancef, Lasix, Mannitol, KCl, crystalloids, colloids, blood products, D5.
*HMC778	Male	65	C	Cerebrovascular Accident	History of alcohol abuse; quit drinking 1972. Possible jail time for drunk and disorderly as youth.	Endocarditis; aortic valve replacement 1987. Rotator cuff surgery 1981, disc surgery 1982, eye surgery 1980. Kidney stones 1998. Hypertension x20 years, controlled by meds. Diabetes x1year, controlled with diet. Mild asthma, no medications. Coumadin, Cardizem, KCl, Anticard. Concussion week prior to admission - kicked in head by horse.	Norcuron, Lasix, Mannitol, Vit K, Dopamine, LCNW, Vasopressin, Insulin, T4, Nipride, blood products, crystalloids
*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.
*HMC786	Male	22	C	Head trauma, MVA	Social alcohol. Positive for marijuana on admission, no social history of drug use.	Biopsy on thyroid nodule 10/01, results negative.	Heparin, Dopamine, crystalloids, tetanus shot, Cefazolin, Clindamycin, Ancef, albumin, Zantac, Mylanta, DDAVP, KCl

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

© DLS

**Donor Information**

Specimen	Gender	Age (years)	Race	Cause of Death	Social History	Medical History	Medication given during Hospitalization
*HMC798	Male	37	C	Cardiopulmonary arrest	Methamphetamines; family states quit 3 months ago, but donor awake 72 hours before being found unresponsive. Jail, 4 months, in past year. Girlfriend also in jail, 2 months, 2001. Social alcohol.	Knee surgery. Elavil.	Heparin, Lasix, Mannitol, Dopamine, Levophed, Epinephrine, Atropine, Bicarb, crystalloids.
HMC800	Male	33	C	Head trauma, ped v. MVA	6 pack beer per day, unknown duration. Marijuana, unknown duration. Multiple tattoos.	Broke right collar bone 2 years ago. Foreign body in right foot, 3 years, never removed. MVA 2 months ago. Allergic to topical ointments. Dog bite 3 years ago; no symptoms.	Heparin, blood products.
*HMC802	Male	61	C	CVA-ICH	Daily beer and whiskey, heavy use. Hospital intake notes alcohol abuse. Nonprofessional tattoos, 1961-62. Jail (days) for DUI, auto theft.	DM, type II, insulin dependent, x10 years, noncompliant with meds. Hypertension x20 years, noncompliant with meds. CAD, history of syncope. MVA in 1969 resulting in head injury.	Heparin, Mannitol, Lidocaine, Dopamine, humulin, accupril, Metoprolol, KCl, insulin, Ancef, Pitressin, Lasix, Solumedrol, blood products.
*HMC810	Male	59	C	CHI s/p fall	4 beers/day. Marijuana (no amount/time provided).	MI 12 years ago, CAD, stent placement 7/02. Coumadin. Cataracts removed 30 years ago. Kidney bilateral, hard aortic plaque.	Heparin, Propofol, Dopamine, Dobutamine, Pentobarb, Vasopressin, Neosynephrne.
HMC819	Male	79	C	CVA	Normal.	Hypertension x5 years, controlled by meds. Carotid study 3 years ago. Cardiac disease, 4 vessels, no treatment. Knee tumor (non cancerous) removed 45 years ago. Glaucoma, resolved spontaneously. Neuro assessment for ALS 1 month ago. Congard, Coumadin, Lozol, vitamins, Selenium, fish oil, ACE with zinc.	Heparin, Dopamine, Kefzol, Vit K, crystalloids.
*HMC829	Male	46	C	Head trauma s/p fall.	4 beers, shot of vodka daily x30 years. Alcohol abuse x10 years. Boyfriend Hep B+ (note: donor's PCR results negative).	Back surgery, bilateral knee surgery s/p MVA, 1997. Weight loss last 6 months. Craniotomy on admission. Allergic to aspirin.	Mannitol, Levophed, Neosynephrine, Solumedrol, DDAVP, insulin, Pavulon, Neomycin, Clindamycin, Nystatin, Erythromycin, Ceftazadime, Ancef, Kphos, KCl, MgSO4, CaGluconate, NaHCO3, crystalloids, colloids, albumin.

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

© DLS



**Donor Information**

Specimen	Gender	Age (years)	Race	Cause of Death	Social History	Medical History	Medication given during Hospitalization
HMC832	Male	72	C	CVA-ICH	Normal.	Hypertension. IDDM x12 years. Valve disease. Mitral valve replacement 1984. Seizures. Schizophrenia. Tardive dyskinesia. Dementia. GI bleed, esophagitis. Allergic rhinitis. Cataract surgery 1995, 2000. Colace, insulin, Coumadin, Lipitor, Depacote, Zantac, Senna, NaCl, Lantus, Kepra, Lisinopril. Allergic to ASA, NSAIDs, Dilantin.	Mannitol, Ativan, Dilantin, Tylenol, Labetolol, Propofol, insulin, Levophed, Dopamine, Kepra, Depakote, Valproic acid, morphine, Unasyn, Lisinopril, Lopressor, Reglan, Nafcillin, Vancomycin, DDAVP, Vecuronium, Etomidate, Succinicholine, Lidocaine, Atracurium, Fentanyl, Norepinephrine, Phenylephrine, KCl, Pepcid, crystalloids.
*HMC834	Male	12	C	Head trauma s/p ATV accident	Normal.	Seizure x1 at 18 months old. Broke collar bone at 3 years old. Vaccinated for Hep B.	Heparin, Mannitol, Levophed, Norepinephrine, Fosphytine, Fentanyl, Versed, Vecuronium, Vancomycin, Cefepime, Pentobarb, TPN, DDAVP, insulin, Pepcid, albumin, blood products.
HMC838	Male	48	C	CVA	12 pack 2x week. Multiple drug use, no injecting, clean 18-20 years. Jail for 3 years and 1 year, 20 years ago, then overnight for DUI 5 years ago.	2 heart attacks in 00. Angiograms in 00 and 03. Lotensin, Plavix, Lopressor, Tricor. Hypertension, controlled with meds. Arthritis, rosacea.	Heparin, Levophed, crystalloids, Decadron, Diprovan, Ms, Ativan, Vasotec, Diltiazem, Levalbutrol, Solumedrol.
*HMC864	Male	69	C	CVA, IPH	Whiskey and beer, a fifth a day for 7-8 years.	Primary emphysema, elevated PSA, hypertension x20 years, benign prostatic hypertrophy, broken ankle. Asbestos exposure, 1960s. BPH. Prostrata, hyperbalance, lung support, herbal meds, digitalis, Coumadin, diocan HC, Saw palmetto.	Heparin, Labetolol, crystalloids, Vecuronium, Versed, Fentanyl, Vitamin K, Mannitol, Lasix
*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.
*HMC911	Male	53	C	CVA 2 <sup>nd</sup> 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.

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

© DLS



**Donor Information**

Specimen	Gender	Age (years)	Race	Cause of Death	Social History	Medical History	Medication given during Hospitalization
HMC912	Male	41	C	Head Trauma-2 <sup>nd</sup> to MVA	Cigarettes;3-4 cigs a day since age of 24. Beer; 2 beers every other day since age 14. Tequila; 2 shots every other day since age 14. Marijuana; Smoked since age 14.	Asthma as a child; outgrown as an adult.	Versed, MGSO4, and Heparin. Flushed and stored in UW solution.
*HMC913	Male	37	C	Anoxia 2 <sup>nd</sup> 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 2 <sup>nd</sup> 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.
HMC928	Male	37	C	ANOXIA 2 <sup>nd</sup> 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 <sup>nd</sup> 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 <sup>nd</sup> 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

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

© DLS

**Donor Information**

Specimen	Gender	Age (years)	Race	Cause of Death	Social History	Medical History	Medication given during Hospitalization
*HMC933	Male	39	C	Head Trauma 2 <sup>nd</sup> 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 2 <sup>nd</sup> 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.

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

© DLS