

H1500.H15T Lot No. 512

Cryopreserved Human Hepatocytes

Donor Information

Demographics

Gender: Female

Age: 62 Years

Race: Caucasian

Cause of Death: Anoxia

Serology

CMV: (-)

HIV: (-)

HBV: (-)

HCV: (-)

Assured Minimum Yield: 4.0×10^6 per vial

Viability: 92.0%

(Yield and viability are based on experiments performed at XenoTech using XenoTech's thawing protocol and K2000 Hepatocyte Isolation Kit.)

Transporter¹

Marker substrate uptake (pmol/million cells/min)^{2§}

| | | |
|------|------------------------------------|-----------|
| OATP | Estrone-3-sulfate | 47 ± 13 |
| NTCP | Taurocholic Acid | 5.5 ± 1.4 |
| OCT1 | 1-Methyl-4-phenylpyridinium Iodine | 18 ± 6 |

Enzyme

Marker substrate reaction (pmol/million cells/min)

| | | |
|----------|--|------|
| CYP2D6 | Dextromethorphan O-demethylation | 66.8 |
| CYP2E1 | Chlorzoxazone 6-hydroxylation | 86.9 |
| CYP3A4/5 | Testosterone 6β-hydroxylation | 43.7 |
| UGT | Glucuronidation of 4-Methylumbelliferone | 366 |

¹ OATP (Organic Anion Transporting Polypeptide), NTCP (Sodium Taurocholate Co-transporting Polypeptide), OCT1 (Organic Cation Transporter)

² Data reflect mean and standard deviation from three separate assays performed on three separate days.

[§] Characterization based on methods described in: Hallifax D and Houston JB (2006) Uptake and intracellular binding of lipophilic amine drugs by isolated rat hepatocytes and implications for prediction of in vivo metabolic clearance. Drug Metabolism and Disposition 34:1829-1836.

CAUTION: These hepatocyte samples are from donors who tested negative for HIV and hepatitis. However, we recommend that these samples be considered as potential biohazards and that universal precautions be used when working with human derived products.

Store vials in liquid nitrogen, vapor phase.



Data sheet prepared 4/21/09