

H1000.H15T **Lot No. H694**

Cryopreserved Human Hepatocytes

Donor Information	
Demographics	Serology
Gender: Male	CMV: (+)
Age: 29 Years	HIV: (-)
Race: Caucasian	HBV: (-)
Cause of Death: Anoxia	HCV: (-)

Assured Minimum Yield: 4.0 x 10⁶ per vial

Viability: 88.6%

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

Transporter ¹	Marker substrate uptake (pmol/million cells/min) ^{2 §}	
OATP	Estrone-3-sulfate	39 ± 12
NTCP	Taurocholic Acid	6.1 ± 1.5
OCT1	1-Methyl-4-phenylpyridinium Iodine	18 ± 5

Enzyme	Marker substrate reaction (pmol/million cells/min)	
CYP2D6	Dextromethorphan O-demethylation	64.6
CYP2E1	Chlorzoxazone 6-hydroxylation	80.0
CYP3A4/5	Testosterone 6β-hydroxylation	14.0
UGT	Glucuronidation of 4-Methylumbelliferone	211

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