

H1500.H15B Lot No. HC4-6

Cryopreserved Human Hepatocytes

Donor Information**Demographics****Gender:** Female**Age:** 56 Years**Race:** Caucasian**Cause of Death:** Cerebrovascular Accident**Serology**

CMV: (+)

HIV: (-)

HTLV: (-)

HBV: (-)

HCV: (-)

Assured Minimum Yield: 4.0×10^6 per vial**Viability:** 78.8%

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

Enzyme**Marker substrate reaction (pmol/million cells/min)**

CYP1A2	Phenacetin <i>O</i> -dealkylation	38.2 ± 5.8
CYP2A6	Coumarin 7-hydroxylation	18.8 ± 1.6
CYP2B6	Bupropion hydroxylation	143 ± 12
CYP2C8	Amodiaquine <i>N</i> -dealkylation	435 ± 31
CYP2C9	Diclofenac 4'-hydroxylation	276 ± 61
CYP2C19	<i>S</i> -Mephenytoin 4'-hydroxylation	2.84 ± 0.49
CYP2D6	Dextromethorphan <i>O</i> -demethylation	60.6 ± 8.9
CYP2E1	Chlorzoxazone 6-hydroxylation	123 ± 18
CYP3A4/5	Testosterone 6 β -hydroxylation	323 ± 60
CYP3A4/5	Midazolam 1'-hydroxylation	39.4 ± 2.4

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 18-Jan-2010