

H1000.H15C Lot No. HC3-3A

Cryopreserved Human Hepatocytes

Donor Information**Demographics**

Gender: Male
Age: 35 Years
Race: Hispanic
Cause of Death: Head Trauma

Serology

CMV: (+)
HIV: (-)
HBV: (-)
HCV: (-)

Assured Minimum Yield: 6.0×10^6 per vial**Viability:** 76.5%

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

Enzyme**Marker substrate reaction (pmol/cells per inc/min)**

CYP1A2	Phenacetin <i>O</i> -dealkylation	53.0 ± 8.0
CYP2A6	Coumarin 7-hydroxylation	13.9 ± 1.1
CYP2B6	Bupropion hydroxylation	17.4 ± 1.9
CYP2C8	Amodiaquine <i>N</i> -dealkylation	114 ± 7
CYP2C9	Diclofenac 4'-hydroxylation	547 ± 58
CYP2C19	<i>S</i> -Mephenytoin 4'-hydroxylation	19.8 ± 2.3
CYP2D6	Dextromethorphan <i>O</i> -demethylation	81.4 ± 3.9
CYP2E1	Chlorzoxazone 6-hydroxylation	238 ± 26
CYP3A4/5	Testosterone 6 β -hydroxylation	83.4 ± 2.1
CYP3A4/5	Midazolam 1'-hydroxylation	14.9 ± 2.5
CYP4A11	Lauric acid 12-hydroxylation	48.9 ± 0.7

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 1/16/09