

H1000.H15B+ Lot No. HC3-15

Cryopreserved Human Hepatocytes

Donor Information

Demographics

Gender: Male
Age: 41 Years
Race: Caucasian
Cause of Death: Anoxia

Serology

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

Assured Minimum Yield: 4.0×10^6 per vial

Viability: 85.0%

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

Enzyme	Marker substrate reaction (pmol/million cells/min)	
CYP1A2	Phenacetin <i>O</i> -dealkylation	7.92 ± 2.93
CYP2A6	Coumarin 7-hydroxylation	16.8 ± 1.4
CYP2B6	Bupropion hydroxylation	6.04 ± 0.14
CYP2C8	Amodiaquine <i>N</i> -dealkylation	100
CYP2C9	Diclofenac 4'-hydroxylation	97.4 ± 5.6
CYP2C19	S-Mephenytoin 4'-hydroxylation	6.95
CYP2D6	Dextromethorphan <i>O</i> -demethylation	5.76
CYP2E1	Chlorzoxazone 6-hydroxylation	BLQ
CYP3A4/5	Testosterone 6β-hydroxylation	31.4
CYP3A4/5	Midazolam 1'-hydroxylation	9.42 ± 1.63
UGT	7-Hydroxycoumarin glucuronidation	553 ± 38
SULT	7-Hydroxycoumarin sulfonation	26.0 ± 1.8

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 10 Feb 2012

H1000.H15B+ Lot No. HC3-15

Cryopreserved Human Hepatocytes

Donor Information

Demographics

Gender: Male
Age: 41 Years
Race: Caucasian
Cause of Death: Anoxia

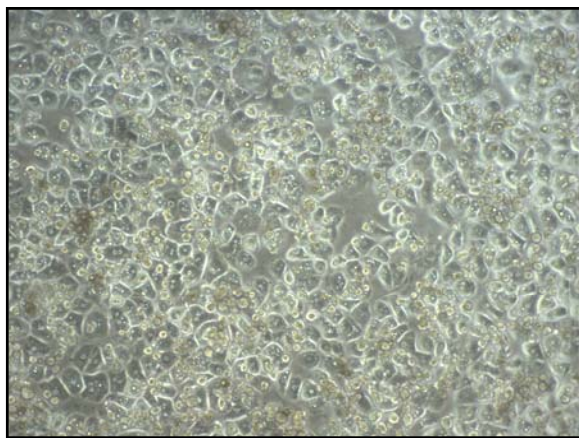
Serology

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

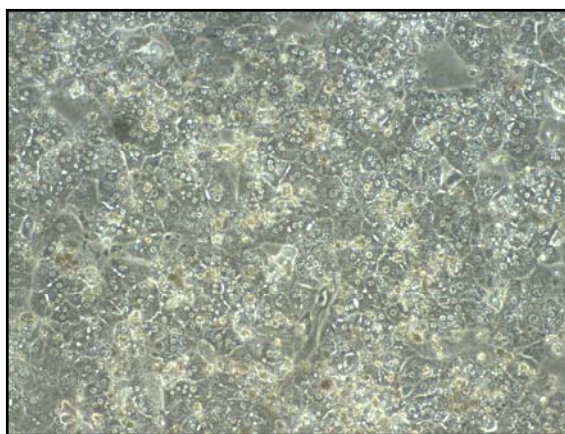
Assured Minimum Yield: 4.0×10^6 per vial

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

Recommended Seeding Density: 1.20×10^6 cells/mL



Photomicrograph (100x) of HC3-15 Day 1 of culture.



Photomicrograph (100x) of HC3-15 prior to incubations.

Induction Data

Enzyme	Prototypical Inducer	Marker substrate reaction	Fold Induction*
CYP1A2	Omeprazole (100 μ M)	Phenacetin O-dealkylation	38.2
CYP2B6	Phenobarbital (750 μ M)	Bupropion hydroxylation	13.6
CYP3A4	Rifampin (10 μ M)	Midazolam 1'-hydroxylation	19.2

* Refers to the increase in pmol metabolite formed when compared to control.

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.

