

H1500.H15C+ Lot No. HC3-51

Cryopreserved Human Hepatocytes
 Human, Female, Individual

Assured Minimum Yield: 6.0×10^6 cells per vial

Viability: 95%

Yield and viability are based on experiments performed at XenoTech using XenoTech's thawing protocol and K8000 OptiThaw Hepatocyte Kit.

Assured Minimum Yield: 6.0×10^6 cells per vial

Viability: 86%

Yield and viability are based on experiments performed at XenoTech using XenoTech's thawing protocol and K2400 (no Percoll gradient) Hepatocyte Isolation Kit.

Enzyme	Marker Substrate Reaction	[S] (μ M)	Rate (pmol/million cells/min)
CYP1A2	Phenacetin O-dealkylation	100	52.5 ± 4.9
CYP2A6	Coumarin 7-hydroxylation	50	62.1 ± 6.5
CYP2B6	Bupropion hydroxylation	500	35.3 ± 1.1
CYP2C8	Amodiaquine N-dealkylation	20	218 ± 6
CYP2C9	Diclofenac 4'-hydroxylation	100	277 ± 34
CYP2C19	S-Mephenytoin 4'-hydroxylation	400	31.9 ± 2.1
CYP2D6	Dextromethorphan O-demethylation	80	44.2 ± 1.5
CYP2E1	Chlorzoxazone 6-hydroxylation	500	121 ± 23
CYP3A4/5	Testosterone 6 β -hydroxylation	250	658 ± 55
CYP3A4/5	Midazolam 1'-hydroxylation	30	78.2 ± 5.0
UGT	7-Hydroxycoumarin glucuronidation	100	605 ± 76
SULT	7-Hydroxycoumarin sulfonation	100	21.2 ± 1.1

Values for enzyme activities were determined at a single substrate concentration and are mean \pm standard deviation of three or more determinations.

To measure cytochrome P450 (CYP), UDP-glucuronosyl transferase (UGT) and sulfotransferase (SULT) activities, hepatocytes (1×10^6 /mL) in suspension were incubated in triplicate at $37 \pm 2^\circ\text{C}$ for 30 minutes in OptiIncubate and marker substrate, at the final concentrations indicated. Metabolite formation was determined by validated LC-MS/MS methods with deuterated metabolites as internal standards.

Donor Information

Gender:	Female
Age:	66 years of age
Race:	Caucasian
Cause of Death:	Head Trauma
Antibody to Cytomegalovirus (CMV):	Negative
All donors tested negative for Human Immunodeficiency Virus (HIV), Hepatitis B Surface Antigen (HBsAg), Hepatitis C Virus, and Rapid Plasma Reagin.	



Store in liquid nitrogen, vapor phase

CAUTION: This sample should be considered as a potential biohazard and universal precautions should be followed. Intended for *in vitro* use only.

These data were generated by and are the property of XenoTech. These data are not to be reproduced, published or distributed without the express written consent of XenoTech.

Datasheet prepared 10 March 2020

H1500.H15C+ Lot No. HC3-51

Cryopreserved Human Hepatocytes
Human, Female, Individual

Assured Minimum Yield: 6.0×10^6 cells per vial

Viability: 95%

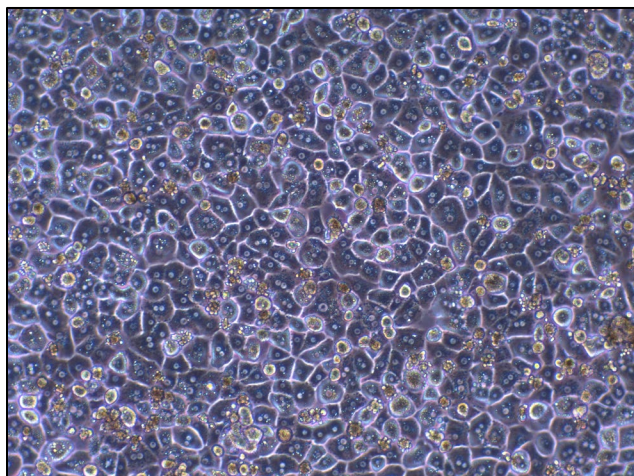
Yield and viability are based on experiments performed at XenoTech using XenoTech's thawing protocol and K8000 OptiThaw Hepatocyte Kit.

Assured Minimum Yield: 6.0×10^6 cells per vial

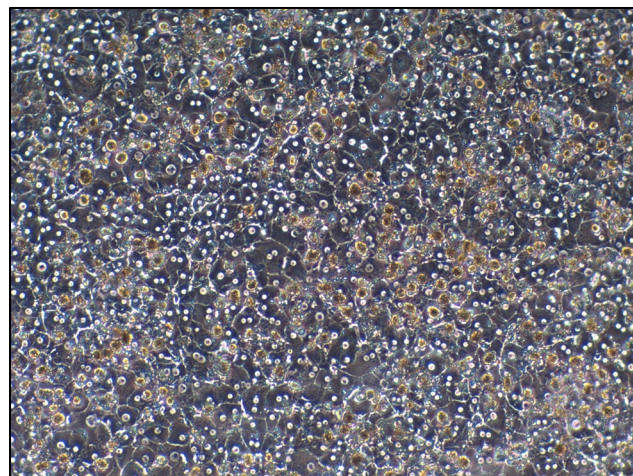
Viability: 86%

Yield and viability are based on experiments performed at XenoTech using XenoTech's thawing protocol and K2400 (no Percoll gradient) Hepatocyte Isolation Kit.

Hepatocyte Cell Culture



Photomicrograph (100x) of HC3-51 Day 1 of culture



Photomicrograph (100x) of HC3-51 incubation day

Plate Format	Recommended Seeding	
	Density (million cells/mL)	Recommended Seeding/ Feeding Volume Per Well
6-well format	1.2	1.7 mL
12-well format	1.2	650 μ L
24-well format	1.2	330 μ L
48-well format	0.75	200 μ L
96-well format	0.75	75 μ L

Induction Data

Enzyme	Inducer	mRNA Fold Induction	Marker Substrate Reaction	Enzymatic Fold Induction
CYP1A2	Omeprazole (50 μ M)	19.6	Phenacetin O-dealkylation	32.5
CYP2B6	Phenobarbital (750 μ M)	4.9	Bupropion hydroxylation	2.5
CYP2B6	CITCO (100 nM)	6.0	Bupropion hydroxylation	3.5
CYP3A4	Rifampin (20 μ M)	5.7	Midazolam 1'-hydroxylation	2.7