

## H1000.H15C+ Lot No. HC0-5

Cryopreserved Human Hepatocytes  
Human, Male, Individual

Assured Minimum Yield:  $6.0 \times 10^6$  cells per vial  
Viability: 86%

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

Enzyme	Marker Substrate Reaction	[S] ( $\mu$ M)	Rate (pmol/million cells/min)
CYP1A2	Phenacetin O-dealkylation	100	177 $\pm$ 10
CYP2A6	Coumarin 7-hydroxylation	50	212 $\pm$ 13
CYP2B6	Bupropion hydroxylation	500	39.0 $\pm$ 3.2
CYP2C8	Amodiaquine N-dealkylation	20	494 $\pm$ 52
CYP2C9	Diclofenac 4'-hydroxylation	100	221 $\pm$ 33
CYP2C19	S-Mephenytoin 4'-hydroxylation	400	28.0 $\pm$ 0.8
CYP2D6	Dextromethorphan O-demethylation	80	22.6 $\pm$ 3.9
CYP2E1	Chlorzoxazone 6-hydroxylation	500	60.5 $\pm$ 2.7
CYP3A4/5	Testosterone 6 $\beta$ -hydroxylation	250	159 $\pm$ 15
CYP3A4/5	Midazolam 1'-hydroxylation	30	154 $\pm$ 18
UGT	7-Hydroxycoumarin glucuronidation	100	560 $\pm$ 11
SULT	7-Hydroxycoumarin sulfonation	100	28.5 $\pm$ 2.7

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 ( $1 \times 10^6$  /mL) in suspension were incubated in triplicate at  $37 \pm 1^\circ\text{C}$  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

<b>Gender:</b>	Male
<b>Age:</b>	78 years of age
<b>Race:</b>	Caucasian
<b>Antibody to Cytomegalovirus (CMV):</b>	Not Determined
<b>Human Immunodeficiency Virus (HIV):</b>	Negative
<b>Hepatitis B Surface Antigen (HbsAg):</b>	Negative
<b>Antibody to Hepatitis C Virus (HCV):</b>	Negative



### 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 03 May 2018

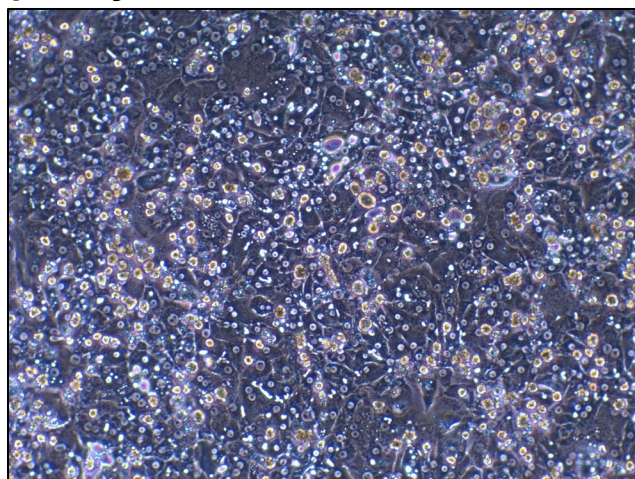
# H1000.H15C+ Lot No. HC0-5

Cryopreserved Human Hepatocytes  
Human, Male, Individual

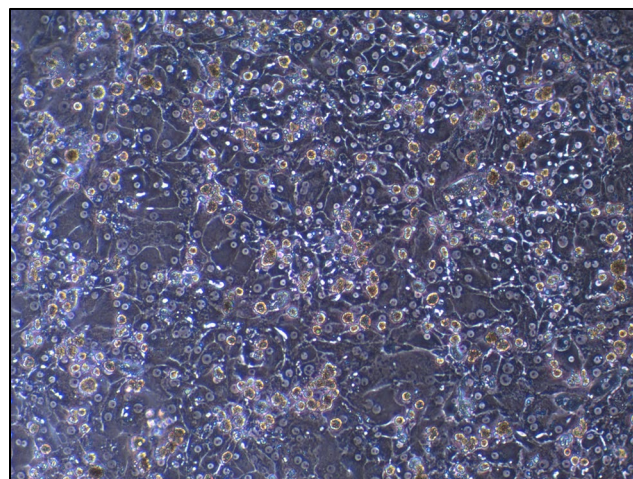
Assured Minimum Yield:  $6.0 \times 10^6$  cells per vial  
Viability: 86%

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

## Hepatocyte Cell Culture



Photomicrograph (100x) of HC0-5 Day 3 of culture



Photomicrograph (100x) of HC0-5 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 $\mu$ L
24-well format	1.2	330 $\mu$ L
48-well format	0.75	200 $\mu$ L
96-well format	0.75	75 $\mu$ L

## Induction Data

Enzyme	Inducer	mRNA Fold Induction	Marker Substrate Reaction	Enzymatic Fold Induction
CYP1A2	Omeprazole (50 $\mu$ M)	240	Phenacetin O-dealkylation	94.2
CYP2B6	Phenobarbital (750 $\mu$ M)	7.2	Bupropion hydroxylation	5.9
CYP2B6	CITCO (100 nM)	8.9	Bupropion hydroxylation	4.7
CYP3A4	Rifampin (20 $\mu$ M)	31.8	Midazolam 1'-hydroxylation	3.7