

## H1000.H15B+ Lot No. HC2-25

Cryopreserved Human Hepatocytes  
Human, Male, Individual

Assured Minimum Yield:  $4.0 \times 10^6$  cells per vial  
Viability: 86.5%

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

Enzyme	Marker Substrate Reaction	[S] ( $\mu$ M)	Rate (pmol/million cells/min)
CYP1A2	Phenacetin O-dealkylation	100	$8.10 \pm 0.48$
CYP2A6	Coumarin 7-hydroxylation	50	$25.3 \pm 7.5$
CYP2B6	Bupropion hydroxylation	500	$105 \pm 18$
CYP2C8	Amodiaquine N-dealkylation	20	$191 \pm 15$
CYP2C9	Diclofenac 4'-hydroxylation	100	$159 \pm 13$
CYP2C19	S-Mephenytoin 4'-hydroxylation	400	$0.667 \pm 0.095$
CYP2D6	Dextromethorphan O-demethylation	80	$26.9 \pm 3.1$
CYP2E1	Chlorzoxazone 6-hydroxylation	500	$167 \pm 8$
CYP3A4/5	Testosterone 6 $\beta$ -hydroxylation	250	$137 \pm 16$
CYP3A4/5	Midazolam 1'-hydroxylation	30	$54.9 \pm 2.1$
UGT	7-Hydroxycoumarin glucuronidation	100	$1350 \pm 210$
SULT	7-Hydroxycoumarin sulfonation	100	$39.8 \pm 6.8$

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 Krebs-Henseleit buffer 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>	58 years of age
<b>Race:</b>	Hispanic
<b>Cause of Death:</b>	Cerebrovascular Accident
<b>Cytomegalovirus (CMV):</b>	Positive
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 24 January 2019

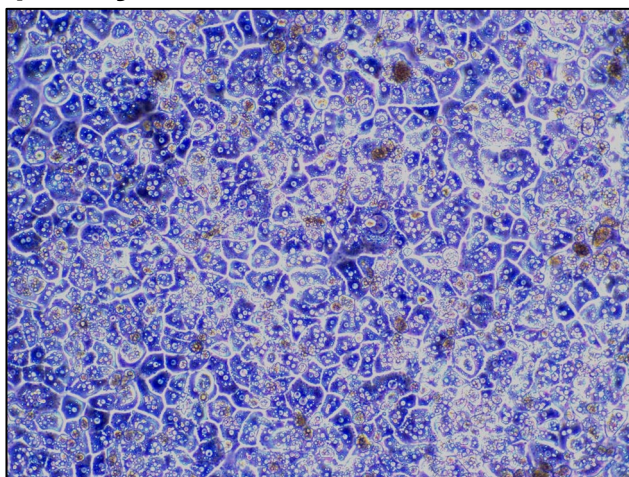
# H1000.H15B+ Lot No. HC2-25

Cryopreserved Human Hepatocytes  
Human, Male, Individual

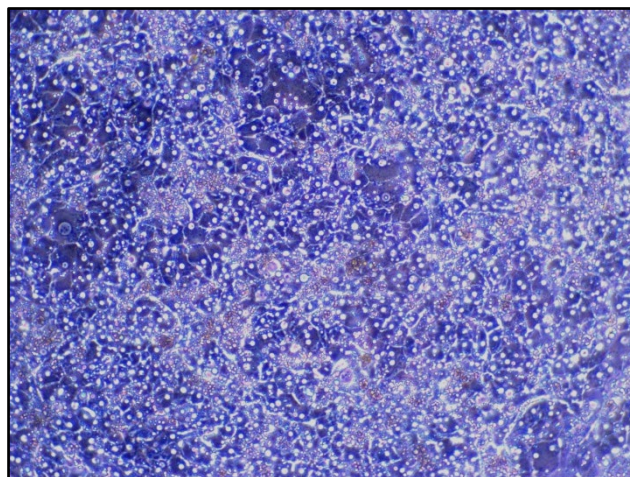
Assured Minimum Yield:  $4.0 \times 10^6$  cells per vial  
Viability: 86.5%

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

## Hepatocyte Cell Culture



Photomicrograph (100x) of HC2-25 Day 2 of culture



Photomicrograph (100x) of HC2-25 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	Not Tested	

## Induction Data

Enzyme	Prototypical Inducer	Marker substrate reaction	Fold Induction*
CYP1A2	Omeprazole (100 $\mu$ M)	Phenacetin O-dealkylation	19.0
CYP2B6	Phenobarbital (750 $\mu$ M)	Bupropion hydroxylation	5.93
CYP3A4	Rifampin (10 $\mu$ M)	Testosterone 6 $\beta$ -hydroxylation	3.71

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