

## H1500.H15C Lot No. HC3-41

Cryopreserved Human Hepatocytes  
Human, Female, Individual

Assured Minimum Yield: 6.0 x 10<sup>6</sup> cells per vial  
Viability: 78%

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

Enzyme	Marker Substrate Reaction	[S] (μM)	Rate (pmol/million cells/min)
CYP1A2	Phenacetin O-dealkylation	100	92.1 ± 12.1
CYP2A6	Coumarin 7-hydroxylation	50	135 ± 20
CYP2B6	Bupropion hydroxylation	500	48.4 ± 2.1
CYP2C8	Amodiaquine N-dealkylation	20	486 ± 20
CYP2C9	Diclofenac 4'-hydroxylation	100	436 ± 33
CYP2C19	S-Mephenytoin 4'-hydroxylation	400	111 ± 9
CYP2D6	Dextromethorphan O-demethylation	80	68.3 ± 5.6
CYP2E1	Chlorzoxazone 6-hydroxylation	500	146 ± 18
CYP3A4/5	Testosterone 6β-hydroxylation	250	1070 ± 170
CYP3A4/5	Midazolam 1'-hydroxylation	30	185 ± 12
UGT	7-Hydroxycoumarin glucuronidation	100	442 ± 58
SULT	7-Hydroxycoumarin sulfonation	100	5.89 ± 0.48

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

To measure cytochrome P450 (CYP), UDP-glucuronosyl transferase (UGT) and sulfotransferase (SULT) activities, hepatocytes (1 x 10<sup>6</sup> /mL) in suspension were incubated in triplicate at 37 ± 1°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>	Female
<b>Age:</b>	56 years of age
<b>Race:</b>	Caucasian
<b>Cause of Death:</b>	Head trauma
<b>Antibody to Cytomegalovirus (CMV):</b>	Positive
<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 14 April 2017