

H1000.H15C Lot No. HC2-47

Cryopreserved Human Hepatocytes
 Human, Male, Individual

Assured Minimum Yield: 6.0×10^6 cells per vial
 Viability: 81%

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

Enzyme	Marker Substrate Reaction	[S] (μ M)	Rate (pmol/million cells/min)
CYP1A2	Phenacetin O-dealkylation	100	76.2 ± 8.0
CYP2A6	Coumarin 7-hydroxylation	50	143 ± 19
CYP2B6	Bupropion hydroxylation	500	127 ± 20
CYP2C8	Amodiaquine N-dealkylation	20	510 ± 8
CYP2C9	Diclofenac 4'-hydroxylation	100	327 ± 44
CYP2C19	S-Mephenytoin 4'-hydroxylation	400	53.4 ± 3.9
CYP2D6	Dextromethorphan O-demethylation	80	69.5 ± 6.2
CYP2E1	Chlorzoxazone 6-hydroxylation	500	45.5 ± 1.8
CYP3A4/5	Testosterone 6 β -hydroxylation	250	643 ± 6
CYP3A4/5	Midazolam 1'-hydroxylation	30	51.2 ± 13.0
UGT	7-Hydroxycoumarin glucuronidation	100	614 ± 32
SULT	7-Hydroxycoumarin sulfonation	100	44.8 ± 2.8

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 (1×10^6 /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

Gender:	Male
Age:	44 years of age
Race:	Caucasian
Cause of Death:	Anoxia
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 09 April 2019