

H1000.H15-B Lot No. HC5-12

Cryopreserved Human Hepatocytes
 Human, Male, Individual

Assured Minimum Yield: 4×10^6 cells per vial
 Average Yield 5.66×10^6 cells per vial
 Average Viability: 77.9%

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] (μM)	Rate ($\mu\text{mol}/\text{million cells}/\text{min}$)
CYP1A2	Phenacetin O-dealkylation	100	19.7 ± 1.5
CYP2A6	Coumarin 7-hydroxylation	50	147 ± 22
CYP2B6	Bupropion hydroxylation	500	199 ± 7
CYP2C8	Amodiaquine N-dealkylation	20	282 ± 32
CYP2C9	Diclofenac 4'-hydroxylation	100	298 ± 72
CYP2C19	S-Mephenytoin 4'-hydroxylation	400	2.19 ± 0.35
CYP2D6	Dextromethorphan O-demethylation	80	48.5 ± 7.6
CYP2E1	Chlorzoxazone 6-hydroxylation	500	239 ± 28
CYP3A4/5	Testosterone 6β -hydroxylation	250	21.7 ± 0.9
CYP3A4/5	Midazolam 1'-hydroxylation	30	16.2 ± 2.0
UGT	7-Hydroxycoumarin glucuronidation	100	462 ± 106
SULT	7-Hydroxycoumarin sulfonation	100	82.9

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 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

Gender:	Male
Age:	56 years
Race:	Caucasian
Cause of Death:	Head trauma
Cytomegalovirus (CMV):	Positive
Human Immunodeficiency Virus (HIV):	Negative
Hepatitis B Surface Antigen (HbsAg):	Negative
Antibody to Hepatitis C Virus (HCV):	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 24 December 2014