

H1000.H15B Lot No. HC3-43

Cryopreserved Human Hepatocytes
Human, Male, Individual

Assured Minimum Yield: 4.0 x 10⁶ cells per vial
Viability: 74%

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	72.4 ± 10.1
CYP2A6	Coumarin 7-hydroxylation	50	35.0 ± 1.9
CYP2B6	Bupropion hydroxylation	500	23.7 ± 3.3
CYP2C8	Amodiaquine N-dealkylation	20	243 ± 22
CYP2C9	Diclofenac 4'-hydroxylation	100	122 ± 4
CYP2C19	S-Mephenytoin 4'-hydroxylation	400	11.0 ± 1.4
CYP2D6	Dextromethorphan O-demethylation	80	14.6 ± 1.2
CYP2E1	Chlorzoxazone 6-hydroxylation	500	195 ± 10
CYP3A4/5	Testosterone 6β-hydroxylation	250	193 ± 13
CYP3A4/5	Midazolam 1'-hydroxylation	30	37.9 ± 2.4
UGT	7-Hydroxycoumarin glucuronidation	100	281 ± 16
SULT	7-Hydroxycoumarin sulfonation	100	10.1 ± 0.5

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 10⁶ /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

Gender:	Male
Age:	61 years of age
Race:	Caucasian
Cause of Death:	Head trauma
Antibody to 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 14 April 2017