

CryostaX

Single Freeze Cryopreserved Human Hepatocytes

HP1500.HP Lot No. H1416

Human, Female, Individual

Assured Minimum Yield: 5.0×10^6 cells per vial
 Viability: 92%

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	92.3 ± 2.6
CYP2A6	Coumarin 7-hydroxylation	50	18.9 ± 2.2
CYP2B6	Bupropion hydroxylation	500	10.8 ± 1.1
CYP2C8	Amodiaquine N-dealkylation	20	140 ± 12
CYP2C9	Diclofenac 4'-hydroxylation	100	68.7 ± 5.3
CYP2C19	S-Mephenytoin 4'-hydroxylation	400	24.5 ± 0.6
CYP2D6	Dextromethorphan O-demethylation	80	20.2 ± 1.6
CYP2E1	Chlorzoxazone 6-hydroxylation	500	70.1 ± 3.4
CYP3A4/5	Testosterone 6 β -hydroxylation	250	290 ± 16
CYP3A4/5	Midazolam 1'-hydroxylation	30	23.8 ± 0.8
UGT	7-Hydroxycoumarin glucuronidation	100	496 ± 31
SULT	7-Hydroxycoumarin sulfonation	100	11.5 ± 0.6

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 $37 \pm 2^\circ\text{C}$ 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:	Female
Age:	67 years of age
Race:	African American
Cause of Death:	Cerebrovascular accident
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 14 January 2021