

CryostaX

Single Freeze Plateable Cryopreserved Human Hepatocytes

HP1000.HP+ Lot No. H1434

Cryopreserved Human Hepatocytes
Human, Male, Individual

Assured Minimum Yield: 5.0×10^6 cells per vial
Viability: 89%

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	42.3 \pm 2.5
CYP2A6	Coumarin 7-hydroxylation	50	45.7 \pm 5.5
CYP2B6	Bupropion hydroxylation	500	41.2 \pm 2.3
CYP2C8	Amodiaquine N-dealkylation	20	456 \pm 45
CYP2C9	Diclofenac 4'-hydroxylation	100	386 \pm 21
CYP2C19	S-Mephenytoin 4'-hydroxylation	400	2.33 \pm 0.27
CYP2D6	Dextromethorphan O-demethylation	80	44.8 \pm 0.8
CYP2E1	Chlorzoxazone 6-hydroxylation	500	96.6 \pm 4.1
CYP3A4/5	Testosterone 6 β -hydroxylation	250	345 \pm 38
CYP3A4/5	Midazolam 1'-hydroxylation	30	105 \pm 11
UGT	7-Hydroxycoumarin glucuronidation	100	882 \pm 45
SULT	7-Hydroxycoumarin sulfonation	100	20.5 \pm 0.1

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:	Male
Age:	55 years of age
Race:	Caucasian
Cause of Death:	Head Trauma
Antibody to Cytomegalovirus (CMV):	Positive
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 19 May 2021

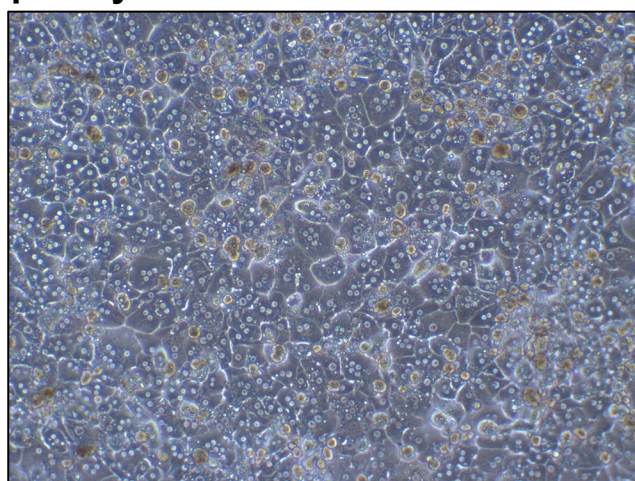
HP1000.HP+ Lot No. H1434

Cryopreserved Human Hepatocytes
Human, Male, Individual

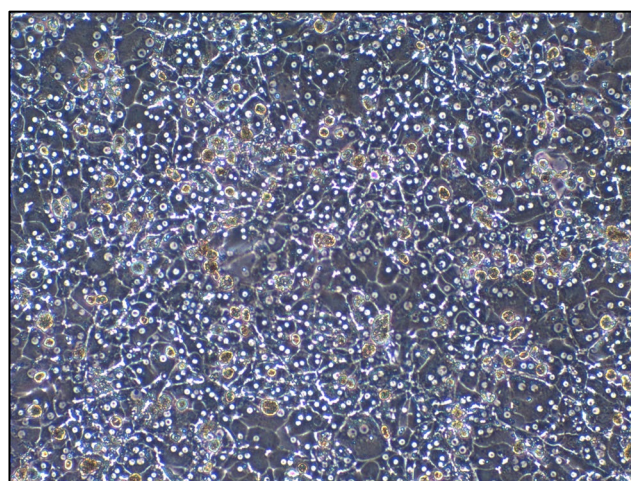
Assured Minimum Yield: 5.0×10^6 cells per vial
Viability: 89%

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

Hepatocyte Cell Culture



Photomicrograph (100x) of H1434 Day 2 of culture



Photomicrograph (100x) of H1434 incubation day

Plate Format	Recommended Seeding	
	Density (million cells/mL)	Recommended Seeding/ Feeding Volume Per Well
6-well format	1.2	1.7 mL
12-well format	1.2	650 μ L
24-well format	1.2	330 μ L
48-well format	0.75	200 μ L
96-well format	0.75	75 μ L

Induction Data

Enzyme	Inducer	mRNA Fold Induction	Marker Substrate Reaction	Enzymatic Fold Induction
CYP1A2	Omeprazole (50 μ M)	172.3	Phenacetin O-dealkylation	100.8
CYP2B6	Phenobarbital (750 μ M)	14.9	Bupropion hydroxylation	4.7
CYP2B6	CITCO (100 nM)	11.5	Bupropion hydroxylation	6.8
CYP3A4	Rifampin (20 μ M)	7.7	Midazolam 1'-hydroxylation	2.1