

CryostaX

Pooled Cryopreserved Human Hepatocytes

HPCH10 Lot No. 2310193

Pool of 10 (5 Female and 5 Male)

Assured Minimum Yield: 5.0 x 10⁶ cells per vial

Viability: 85%

Enzyme	Marker Substrate Reaction	[S] (µM)	Rate (pmol/million cells/min)
CYP1A2	Phenacetin O-dealkylation	100	95.1 ± 1.1
CYP2A6	Coumarin 7-hydroxylation	50	28.5 ± 3.6
CYP2B6	Bupropion hydroxylation	500	33.2 ± 2.6
CYP2C8	Amodiaquine N-dealkylation	20	346 ± 51
CYP2C9	Diclofenac 4'-hydroxylation	100	241 ± 3
CYP2C19	S-Mephenytoin 4'-hydroxylation	400	40.3 ± 1.1
CYP2D6	Dextromethorphan O-demethylation	80	47.5 ± 1.7
CYP2E1	Chlorzoxazone 6-hydroxylation	500	128 ± 16
CYP3A4/5	Testosterone 6β-hydroxylation	250	402 ± 8
CYP3A4/5	Midazolam 1'-hydroxylation	30	113 ± 15
UGT	7-Hydroxycoumarin glucuronidation	100	676 ± 11
SULT	7-Hydroxycoumarin sulfonation	100	20.3 ± 0.8

To measure cytochrome P450 (CYP), UDP-glucuronosyl transferase (UGT) and sulfotransferase (SULT) activities, hepatocytes (1 x 10^6 cells/mL) in suspension were incubated in triplicate at $37 \pm 2^{\circ}$ C for 30 minutes in Optilncubate and marker substrate, at the final concentrations indicated. Metabolite formation was determined by validated LC-MS/MS methods with deuterated metabolites as internal standards.

Uptake Activity Data

Uptake Transporter	Marker Substrate	[S] (µM)	Rate (pmol/million cells/min)
OATP1B1	Estrone sulfate	1	18.7
OATP1B3	CCK-8	1	6.1
OCT1	MPP+	1	6.0
NTCP	TCA	1	4.2

To measure uptake activities, hepatocytes (1.0 x 10⁶ cells/mL) in suspension were incubated in triplicate at 4°C and 37°C for 1 minute in Krebs-Henseleit buffer and marker substrate, at the final concentrations indicated. Uptake of substrate was measured by scintillation counter.

Donor Information

Gender: Males (5), Females (5) **Age:** 33-67 years of age

Race: African American (2), Caucasian (8)

Cause of Death: Anoxia (2), Cerebrovascular accident (5), Head trauma (3)

Antibody to Cytomegalovirus (CMV): Positive (6), Negative (4)

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.

This data sheet serves as a Certificate of Analysis and has been approved by Stephanie Helmstetter, Assistant Director.

Signature and Date:

Assistant Director.

15 August 2023