

# CryostaX

Single-Freeze Pooled Cryopreserved Human Hepatocytes

**HPCH20-50**

**Lot No. 2310103**

Pool of 20

Assured Minimum Yield:

5.0 x 10<sup>6</sup> cells per vial

Viability:

74%

Enzyme	Marker Substrate Reaction	[S] (μM)	Rate (pmol/million cells/min)
CYP1A2	Phenacetin O-dealkylation	100	71.9 ± 4.1
CYP2A6	Coumarin 7-hydroxylation	50	67.3 ± 1.0
CYP2B6	Bupropion hydroxylation	500	84.8 ± 5.7
CYP2C8	Amodiaquine N-dealkylation	20	476 ± 36
CYP2C9	Diclofenac 4'-hydroxylation	100	264 ± 12
CYP2C19	S-Mephenytoin 4'-hydroxylation	400	10.4 ± 0.6
CYP2D6	Dextromethorphan O-demethylation	80	65.1 ± 0.5
CYP2E1	Chlorzoxazone 6-hydroxylation	500	141 ± 3
CYP3A4/5	Testosterone 6β-hydroxylation	250	333 ± 15
CYP3A4/5	Midazolam 1'-hydroxylation	30	67.5 ± 3.0
UGT	7-Hydroxycoumarin glucuronidation	100	478 ± 25
SULT	7-Hydroxycoumarin sulfonation	100	22.7 ± 1.4

To measure cytochrome P450 (CYP), UDP-glucuronosyl transferase (UGT) and sulfotransferase (SULT) activities, hepatocytes (1 x 10<sup>6</sup> cells/mL) in suspension were incubated in triplicate at 37 ± 2°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.

## Uptake Activity Data

Uptake Transporter	Marker Substrate	[S] (μM)	Rate (pmol/million cells/min)
OATP1B1	Estrone sulfate	1	27.0
OATP1B3	CCK-8	1	7.6
OCT1	MPP+	1	3.1
NTCP	TCA	1	4.8

To measure uptake activities, hepatocytes (0.5 x 10<sup>6</sup> cells/mL) in suspension were incubated in triplicate at 4°C ± 2°C and 37°C ± 2°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

<b>Gender:</b>	Males (10), Females (10)
<b>Age:</b>	7-65 years of age
<b>Race:</b>	Caucasian (18), African American (2)
<b>Cause of Death:</b>	Cerebrovascular accident (7), Anoxia (11), Head trauma (2)
<b>Antibody to Cytomegalovirus (CMV):</b>	Positive (10), Negative (9), Not Determined (1)
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: Stephanie Helmstetter 19 April 2023