

# CryostaX Geneknown™

Single Freeze Pooled Cryopreserved Human Hepatocytes

## HPCH.UGT1A1.HA Lot No. 1510253

UGT1A1 Genotype: \*1/\*1

Pool of 3

Assured Minimum Yield:  $4.5 \times 10^6$  cells per vial

Viability: >70.0%

### Individual Donor CYP Genotype Information:

Donor	CYP1A1	CYP1A2	CYP2A6	CYP2B6	CYP2D6	CYP2C8	CYP2C9	CYP2C19	CYP2E1	CYP3A4	CYP3A5
1022	*1/*1	*1/*1F	*1/*1	*1/*8	*3/*41	*1/*4	*1/*1	*1/*1	*1/*1	*1/*1	*3/*3
1197	*1/*1	*1/*1F	*1/*1	*1/*6	*1/*2	*1/*1	*1/*1	*1/*1	*1/*1	*1/*22	*3/*3
1208	*1/*1	*1/*1	*1/*1	*1/*1	*1/*1	*1/*1	*1/*3	*1/*1	*1/*1	*1/*1	*3/*3

Enzyme	Marker Substrate Reaction	[S] (μM)	Rate (pmol/million cells/min)
CYP1A2	Phenacetin O-dealkylation	100	62.9
CYP2A6	Coumarin 7-hydroxylation	50	31.4
CYP2B6	Bupropion hydroxylation	500	19.2
CYP2C8	Amodiaquine N-dealkylation	20	88.1
CYP2C9	Diclofenac 4'-hydroxylation	100	209
CYP2C19	S-Mephenytoin 4'-hydroxylation	400	15.4
CYP2D6	Dextromethorphan O-demethylation	80	26.1
CYP2E1	Chlorzoxazone 6-hydroxylation	500	47.4
CYP3A4/5	Testosterone 6β-hydroxylation	250	130
CYP3A4/5	Midazolam 1'-hydroxylation	30	31.4

Values for enzyme activities are averages of the individual donor activities for each specified CYP, therefore the values indicated should be considered theoretical.

To measure cytochrome P450 (CYP) activities, hepatocytes ( $1 \times 10^6$  cells/mL) in suspension were incubated in triplicate at  $37 \pm 1^\circ\text{C}$  for 30 minutes in Krebs-Henseleit buffer 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

<b>Gender:</b>	Male (1), Female (2)
<b>Age:</b>	41-67 years of age
<b>Race:</b>	Caucasian (3)
<b>Cause of Death:</b>	Anoxia (2), Cerebrovascular accident (1)
<b>Cytomegalovirus (CMV):</b>	Positive (3)
<b>Human Immunodeficiency Virus (HIV):</b>	Negative (3)
<b>Hepatitis B Surface Antigen (HBsAg):</b>	Negative (3)
<b>Antibody to Hepatitis C Virus (HCV):</b>	Negative (3)



### 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 26 January 2016