

H1500.H15C+ Lot No. HC5-29

Cryopreserved Human Hepatocytes
 Human, Female, Individual

Assured Minimum Yield: 6.0×10^6 cells per vial
 Viability: 85.0%

Yield and viability are based on experiments performed at XenoTech using XenoTech's thawing protocol and K2000 Hepatocyte Isolation Kit.

Enzyme	Marker Substrate Reaction	[S] (μM)	Rate (pmol/million cells/min)
CYP1A2	Phenacetin O-dealkylation	100	7.69 ± 0.23
CYP2A6	Coumarin 7-hydroxylation	50	24.1 ± 4.0
CYP2B6	Bupropion hydroxylation	500	21.5 ± 0.5
CYP2C8	Amodiaquine N-dealkylation	20	171 ± 17
CYP2C9	Diclofenac 4'-hydroxylation	100	199 ± 32
CYP2C19	S-Mephenytoin 4'-hydroxylation	400	2.50 ± 0.27
CYP2D6	Dextromethorphan O-demethylation	80	23.5 ± 1.9
CYP2E1	Chlorzoxazone 6-hydroxylation	500	197 ± 12
CYP3A4/5	Testosterone 6β -hydroxylation	250	94.2 ± 18.2
CYP3A4/5	Midazolam 1'-hydroxylation	30	47.6 ± 10.0
UGT	7-Hydroxycoumarin glucuronidation	100	584 ± 6
SULT	7-Hydroxycoumarin sulfonation	100	26.9 ± 0.9

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 $37 \pm 1^\circ\text{C}$ 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

Gender:	Female
Age:	69 years of age
Race:	Asian
Cause of Death:	Cerebrovascular Accident
Cytomegalovirus (CMV):	Positive
Human Immunodeficiency Virus (HIV):	Negative
Hepatitis B Surface Antigen (HbsAg):	Negative
Antibody to Hepatitis C Virus (HCV):	Negative



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 16 July 2014

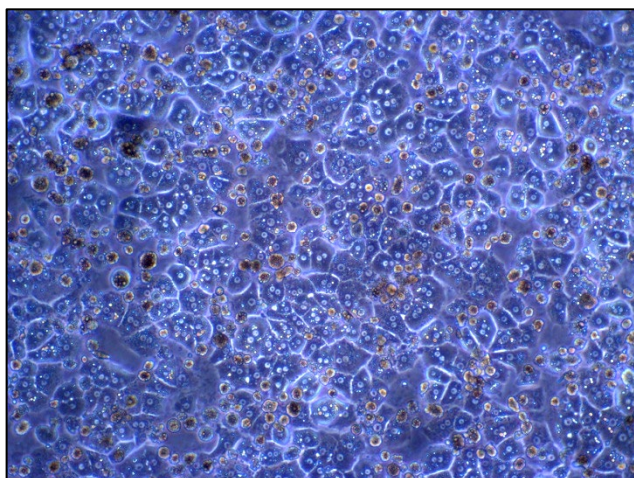
H1500.H15C+ Lot No. HC5-29

Cryopreserved Human Hepatocytes
Human, Female, Individual

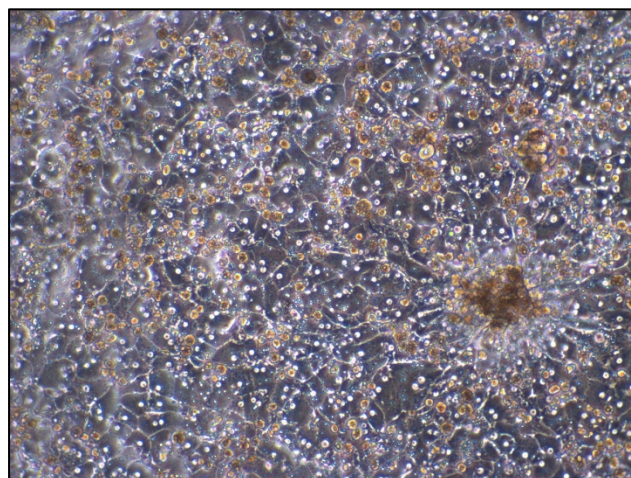
Assured Minimum Yield: 6.0×10^6 cells per vial
Viability: 85.0%

Yield and viability are based on experiments performed at XenoTech using XenoTech's thawing protocol and K2000 Hepatocyte Isolation Kit.

Hepatocyte Cell Culture



Photomicrograph (100x) of HC5-29 Day 2 of culture



Photomicrograph (100x) of HC5-29 incubation day

Plate Format	Recommended Seeding	
	Density (million cells/mL)	Recommended Seeding/ Feeding Volume Per Well
6-well format	1.4	1.7 mL
12-well format	1.4	650 μ L
24-well format	1.4	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 (100 μ M)	30.3	Phenacetin O-dealkylation	47.2
CYP2B6	Phenobarbital (750 μ M)	8.54	Bupropion hydroxylation	19.1
CYP2B6	CITCO (100 nM)	4.88	Bupropion hydroxylation	Not determined
CYP3A4	Rifampin (10 μ M)	14.6	Testosterone 6 β -hydroxylation	4.37