

HU1A1.HA Lot No. 0710412

Human Liver Microsomes

Male, Individual No. 509

0.5 mL at 20 mg protein / mL

Genotype, specific content and activities ^a Content / Rate

UGT1A1 Allelic variant	UGT1A1*1/*1
Cytochrome P450 (nmol/mg protein)	0.243
Cytochrome b ₅ (nmol/mg protein)	0.344
NADPH-cytochrome c reductase (nmol/mg protein/min)	153 ± 3

Enzyme	Marker substrate reaction (pmol/mg protein/min)	
CYP1A2	Phenacetin <i>O</i> -dealkylation	545 ± 23
CYP2A6	Coumarin 7-hydroxylation	272 ± 10
CYP2B6	Bupropion hydroxylation	102 ± 4
CYP2C8	Amodiaquine <i>N</i> -dealkylation	470 ± 8
CYP2C9	Diclofenac 4'-hydroxylation	1690 ± 50
CYP2C19	<i>S</i> -Mephenytoin 4'-hydroxylation	150 ± 5
CYP2D6	Dextromethorphan <i>O</i> -demethylation	155 ± 4
CYP2E1	Chlorzoxazone 6-hydroxylation	2990 ± 180
CYP3A4/5	Testosterone 6β-hydroxylation	571 ± 48
CYP3A4	Midazolam 1'-hydroxylation	36.4 ± 2.5
CYP4A11	Lauric acid 12-hydroxylation	1390 ± 40
UGT1A1	17β-Estradiol 3-glucuronidation	623 ± 4

^a Values for enzyme activities are mean ± standard deviation of three or more determinations.

Sample	Gender	Age (yrs)	Race	Cause of Death
H0509	Male	32	Caucasian	Head Trauma

Serology information

- This donor tested negative for cytomegalovirus
- This donor tested negative for HIV, HTLV, HbsAg, and HCV*
- This donor tested negative for RPR**

* Antibody to Human Immunodeficiency Virus, Antibody to Human T Cell Lymphotropic Virus, Hepatitis B Surface Antigen, Antibody to Hepatitis C Virus, respectively.

** Rapid Plasma Reagin.



Store at -80°C

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, LLC. These data are not to be reproduced, published or distributed without the express written consent of XENOTECH, LLC

DATASHEET PREPARED 20-APRIL-10