

HU1A1.NA Lot No. 0710411

Human Liver Microsomes

Female, Individual No. 508

0.5 mL at 20 mg protein / mL

Genotype, specific content and activities ^a		Content / Rate
UGT1A1 Allelic variant		UGT1A1*28/*28
Cytochrome P450 (nmol/mg protein)		0.336
Cytochrome b ₅ (nmol/mg protein)		0.254
NADPH-cytochrome c reductase (nmol/mg protein/min)		98.9 ± 0.7
Enzyme	Marker substrate reaction (pmol/mg protein/min)	
CYP1A2	Phenacetin <i>O</i> -dealkylation	599 ± 28
CYP2A6	Coumarin 7-hydroxylation	80.4 ± 5.7
CYP2B6	Bupropion hydroxylation	124 ± 5
CYP2C8	Amodiaquine <i>N</i> -dealkylation	1120 ± 40
CYP2C9	Diclofenac 4'-hydroxylation	1270 ± 20
CYP2C19	<i>S</i> -Mephenytoin 4'-hydroxylation	21.5 ± 2.0
CYP2D6	Dextromethorphan <i>O</i> -demethylation	16.6 ± 1.1
CYP2E1	Chlorzoxazone 6-hydroxylation	1410 ± 180
CYP3A4/5	Testosterone 6β-hydroxylation	2250 ± 340
CYP3A4	Midazolam 1'-hydroxylation	53.9 ± 1.8
CYP4A11	Lauric acid 12-hydroxylation	1460 ± 70
UGT1A1	17β-Estradiol 3-glucuronidation	210 ± 3

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

Sample	Gender	Age (yrs)	Race	Cause of Death
H0508	Female	58	Caucasian	Cerebrovascular Accident

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