

HU1A9.HA Lot No. 0910052

Human Liver Microsomes

Male, Individual No. 440

0.5 mL at 20 mg protein / mL

Suspension medium: 250 mM sucrose

Genotype, specific content and activities ^a Content / Rate

UGT1A9 Allelic variant	UGT1A9*1/*1
Cytochrome P450 (nmol/mg protein)	0.118
Cytochrome b ₅ (nmol/mg protein)	0.204
NADPH-cytochrome c reductase (nmol/mg protein/min)	164 ± 0

Enzyme	Marker substrate reaction (pmol/mg protein/min)	Content / Rate
CYP1A2	Phenacetin O-dealkylation	51.5 ± 7.0
CYP2A6	Coumarin 7-hydroxylation	16.3 ± 0.5
CYP2B6	Bupropion hydroxylation	18.5 ± 2.7
CYP2C8	Amodiaquine N-dealkylation	29.7 ± 1.3
CYP2C9	Diclofenac 4'-hydroxylation	844 ± 12
CYP2C19	S-Mephenytoin 4'-hydroxylation	5.25 ± 0.10
CYP2D6	Dextromethorphan O-demethylation	148 ± 5
CYP2E1	Chlorzoxazone 6-hydroxylation	1380 ± 70
CYP3A4/5	Testosterone 6β-hydroxylation	80.7 ± 6.1
CYP3A4	Midazolam 1'-hydroxylation	28.2 ± 0.6
CYP4A11	Lauric acid 12-hydroxylation	524 ± 9
UGT1A9	Propofol glucuronidation	1110 ± 30

^a Values for enzyme activities were determined at a single substrate concentration and are mean ± standard deviation of three or more determinations.

Sample	Gender	Age (yrs)	Race	Cause of Death
H0440	Male	67	Caucasian	Anoxia

Serology information

- This donor tested positive 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 09-JULY-09