

H2C9.MA / Lot No. 0710427

Human Liver Microsomes

Female, Individual No. 538

0.5 mL at 20 mg protein / mL

Genotype, specific content and activities ^a		Content / Rate
CYP2C9 Allelic variant		CYP2C9*1/*2
Cytochrome P450 (nmol/mg protein)		0.470
Cytochrome b ₅ (nmol/mg protein)		0.349
NADPH-cytochrome c reductase (nmol/mg protein/min)		159 ± 6
Enzyme	Marker substrate reaction (pmol/mg protein/min)	
CYP1A2	Phenacetin O-dealkylation	454 ± 47
CYP2A6	Coumarin 7-hydroxylation	1640 ± 90
CYP2B6	Bupropion hydroxylation	284 ± 5
CYP2C8	Amodiaquine N-dealkylation	1060 ± 60
CYP2C9	Diclofenac 4'-hydroxylation	3650 ± 260
CYP2C19	S-Mephenytoin 4'-hydroxylation	72.8 ± 4.1
CYP2D6	Dextromethorphan O-demethylation	141 ± 6
CYP2E1	Chlorzoxazone 6-hydroxylation	2140 ± 110
CYP3A4/5	Testosterone 6β-hydroxylation	3780 ± 480
CYP3A4	Midazolam 1'-hydroxylation	321 ± 14
CYP4A11	Lauric acid 12-hydroxylation	1950 ± 30

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

Sample	Gender	Age (yrs)	Race	Cause of Death
H0538	Female	78	Caucasian	Cerebrovascular accident

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.

Data sheet prepared 2/28/08



Store at -80 °C

For in vitro use only

CAUTION: This liver sample is from a donor who tested negative for HIV and hepatitis. However, we recommend that these samples be considered as potential biohazards and that universal precautions be used when working with human derived products.

These data were generated by and are the property of XENOTECH, LLC. These data are not to be reproduced, published or distributed without the expressed written consent of XENOTECH, LLC.