

H2D6.MA / Lot No. 0710415

Human Liver Microsomes

Female, Individual No. 512

0.5 mL at 20 mg protein / mL

Genotype, specific content and activities ^a		Content / Rate
CYP2D6 Allelic variant		CYP2D6*2/*2
Cytochrome P450 (nmol/mg protein)		0.193
Cytochrome b ₅ (nmol/mg protein)		0.254
NADPH-cytochrome c reductase (nmol/mg protein/min)		111 ± 4
Enzyme	Marker substrate reaction (pmol/mg protein/min)	
CYP1A2	Phenacetin O-dealkylation	294 ± 34
CYP2A6	Coumarin 7-hydroxylation	125 ± 3
CYP2B6	Bupropion hydroxylation	104 ± 6
CYP2C8	Amodiaquine N-dealkylation	268 ± 3
CYP2C9	Diclofenac 4'-hydroxylation	1110 ± 110
CYP2C19	S-Mephenytoin 4'-hydroxylation	6.64 ± 0.41
CYP2D6	Dextromethorphan O-demethylation	191 ± 2
CYP2E1	Chlorzoxazone 6-hydroxylation	686 ± 30
CYP3A4/5	Testosterone 6β-hydroxylation	1230 ± 80
CYP3A4	Midazolam 1'-hydroxylation	67.7 ± 4.6
CYP4A11	Lauric acid 12-hydroxylation	740 ± 39

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

Sample	Gender	Age (yrs)	Race	Cause of Death
H0512	Female	62	Caucasian	Anoxia

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.

Data sheet prepared 2/12/08



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.