

H2D6.MA / Lot No. 0710455

Human Liver Microsomes

Female, Individual No. 443

0.5 mL at 20 mg protein / mL

Genotype, specific content and activities ^a		Content / Rate
CYP2D6 Allelic variant		CYP2D6*1/*4
Cytochrome P450 (nmol/mg protein)		0.327
Cytochrome b ₅ (nmol/mg protein)		0.229
NADPH-cytochrome c reductase (nmol/mg protein/min)		89.8 ± 6.1
Enzyme	Marker substrate reaction (pmol/mg protein/min)	
CYP1A2	Phenacetin O-dealkylation	106 ± 17
CYP2A6	Coumarin 7-hydroxylation	132 ± 12
CYP2B6	Bupropion hydroxylation	73.0 ± 5.2
CYP2C8	Amodiaquine N-dealkylation	256 ± 8
CYP2C9	Diclofenac 4'-hydroxylation	1410 ± 70
CYP2C19	S-Mephenytoin 4'-hydroxylation	5.17 ± 0.47
CYP2D6	Dextromethorphan O-demethylation	98.6 ± 5.7
CYP2E1	Chlorzoxazone 6-hydroxylation	941 ± 21
CYP3A4/5	Testosterone 6β-hydroxylation	867 ± 159
CYP3A4	Midazolam 1'-hydroxylation	173 ± 8
CYP4A11	Lauric acid 12-hydroxylation	624 ± 31

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

Sample	Gender	Age (yrs)	Race	Cause of Death
H0443	Female	64	Hispanic	Cerebrovascular Stroke

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 2/6/07



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.