

**H2D6.MA / Lot No. 0710450**

Human Liver Microsomes

Male, Individual No. 397

0.5 mL at 20 mg protein / mL

**Genotype, specific content and activities <sup>a</sup> Content / Rate**

CYP2D6 Allelic variant	CYP2D6*2/*41
Cytochrome P450 (nmol/mg protein)	0.629
Cytochrome b <sub>5</sub> (nmol/mg protein)	0.364
NADPH-cytochrome c reductase (nmol/mg protein/min)	169 ± 10

Enzyme	Marker substrate reaction (pmol/mg protein/min)	
CYP1A2	Phenacetin <i>O</i> -dealkylation	638 ± 15
CYP2A6	Coumarin 7-hydroxylation	2720 ± 50
CYP2B6	Bupropion hydroxylation	3200 ± 140
CYP2C8	Amodiaquine <i>N</i> -dealkylation	4710 ± 100
CYP2C9	Diclofenac 4'-hydroxylation	2880 ± 60
CYP2C19	<i>S</i> -Mephenytoin 4'-hydroxylation	81.3 ± 2.3
CYP2D6	Dextromethorphan <i>O</i> -demethylation	312 ± 3
CYP2E1	Chlorzoxazone 6-hydroxylation	1790 ± 30
CYP3A4/5	Testosterone 6β-hydroxylation	7680 ± 30
CYP3A4	Midazolam 1'-hydroxylation	1010 ± 80
CYP4A11	Lauric acid 12-hydroxylation	1400 ± 50

<sup>a</sup> Values for enzyme activities are mean ± standard deviation of three or more determinations.

Sample	Gender	Age (yrs)	Race	Cause of Death
H0397	Male	60	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.

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.