

## H2D6.MA / Lot No. 0710446

Human Liver Microsomes

Female, Individual No. 500

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)	0.285
Cytochrome b <sub>5</sub> (nmol/mg protein)	0.319
NADPH-cytochrome c reductase (nmol/mg protein/min)	149 ± 2

Enzyme	Marker substrate reaction (pmol/mg protein/min)	
CYP1A2	Phenacetin <i>O</i> -dealkylation	147 ± 5
CYP2A6	Coumarin 7-hydroxylation	681 ± 27
CYP2B6	Bupropion hydroxylation	77.6 ± 5.6
CYP2C8	Amodiaquine <i>N</i> -dealkylation	658 ± 50
CYP2C9	Diclofenac 4'-hydroxylation	795 ± 115
CYP2C19	<i>S</i> -Mephenytoin 4'-hydroxylation	28.9 ± 2.6
CYP2D6	Dextromethorphan <i>O</i> -demethylation	138 ± 3
CYP2E1	Chlorzoxazone 6-hydroxylation	2970 ± 170
CYP3A4/5	Testosterone 6β-hydroxylation	585 ± 23
CYP3A4	Midazolam 1'-hydroxylation	20.5 ± 0.8
CYP4A11	Lauric acid 12-hydroxylation	980 ± 25

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

Sample	Gender	Age (yrs)	Race	Cause of Death
H0500	Female	51	Caucasian	Head trauma

#### 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 2/27/08



## 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.