

## H2D6.MA / Lot No. 0710457

Human Liver Microsomes  
 Female, Individual No. 454  
 0.5 mL at 20 mg protein / mL

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

CYP2D6 Allelic variant	CYP2D6*1/*3
Cytochrome P450 (nmol/mg protein)	0.360
Cytochrome b <sub>5</sub> (nmol/mg protein)	0.244
NADPH-cytochrome c reductase (nmol/mg protein/min)	117 ± 6

Enzyme	Marker substrate reaction (pmol/mg protein/min)	
CYP1A2	Phenacetin <i>O</i> -dealkylation	119 ± 18
CYP2A6	Coumarin 7-hydroxylation	1390 ± 110
CYP2B6	Bupropion hydroxylation	943 ± 44
CYP2C8	Amodiaquine <i>N</i> -dealkylation	1500 ± 30
CYP2C9	Diclofenac 4'-hydroxylation	1410 ± 50
CYP2C19	<i>S</i> -Mephenytoin 4'-hydroxylation	33.2 ± 1.5
CYP2D6	Dextromethorphan <i>O</i> -demethylation	188 ± 3
CYP2E1	Chlorzoxazone 6-hydroxylation	582 ± 70
CYP3A4/5	Testosterone 6β-hydroxylation	3150 ± 240
CYP3A4	Midazolam 1'-hydroxylation	244 ± 16
CYP4A11	Lauric acid 12-hydroxylation	1140 ± 40

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

Sample	Gender	Age (yrs)	Race	Cause of Death
H0454	Female	59	Caucasian	Cerebrovascular accident

#### 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 2/28/08



## 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.