

## H2D6.NA / Lot No. 0810010

Human Liver Microsomes

Male, Individual No. 502

0.5 mL at 20 mg protein / mL

Genotype, specific content and activities <sup>a</sup>		Content / Rate
CYP2D6 Allelic variant		CYP2D6*4/*4
Cytochrome P450 (nmol/mg protein)		0.160
Cytochrome b <sub>5</sub> (nmol/mg protein)		0.184
NADPH-cytochrome c reductase (nmol/mg protein/min)		102 ± 1
Enzyme	Marker substrate reaction (pmol/mg protein/min)	
CYP1A2	Phenacetin O-dealkylation	240 ± 24
CYP2A6	Coumarin 7-hydroxylation	431 ± 8
CYP2B6	Bupropion hydroxylation	98.7 ± 2.0
CYP2C8	Amodiaquine N-dealkylation	569 ± 29
CYP2C9	Diclofenac 4'-hydroxylation	495 ± 21
CYP2C19	S-Mephenytoin 4'-hydroxylation	1.24 ± 0.15
CYP2D6	Dextromethorphan O-demethylation	10.8 ± 0.9
CYP2E1	Chlorzoxazone 6-hydroxylation	1090 ± 90
CYP3A4/5	Testosterone 6β-hydroxylation	572 ± 6
CYP3A4	Midazolam 1'-hydroxylation	19.9 ± 1.0
CYP4A11	Lauric acid 12-hydroxylation	754 ± 15

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

Sample	Gender	Age (yrs)	Race	Cause of Death
H0502	Male	49	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 2/12/08



## 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.