

## H2D6.NA / Lot No. 0710443

Human Liver Microsomes

Female, Individual No. 486

0.5 mL at 20 mg protein / mL

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

### Content / Rate

CYP2D6 Allelic variant

CYP2D6\*4/\*5

Cytochrome P450 (nmol/mg protein)

0.503

Cytochrome b<sub>5</sub> (nmol/mg protein)

0.314

NADPH-cytochrome c reductase (nmol/mg protein/min)

172 ± 2

#### Enzyme

#### Marker substrate reaction (pmol/mg protein/min)

CYP1A2	Phenacetin <i>O</i> -dealkylation	357 ± 17
CYP2A6	Coumarin 7-hydroxylation	1550 ± 50
CYP2B6	Bupropion hydroxylation	2240 ± 70
CYP2C8	Amodiaquine <i>N</i> -dealkylation	2520 ± 70
CYP2C9	Diclofenac 4'-hydroxylation	3620 ± 50
CYP2C19	<i>S</i> -Mephenytoin 4'-hydroxylation	29.4 ± 2.4
CYP2D6	Dextromethorphan <i>O</i> -demethylation	55.2 ± 1.3
CYP2E1	Chlorzoxazone 6-hydroxylation	3900
CYP3A4/5	Testosterone 6β-hydroxylation	2290 ± 60
CYP3A4	Midazolam 1'-hydroxylation	183 ± 10
CYP4A11	Lauric acid 12-hydroxylation	2070 ± 60

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

Sample	Gender	Age (yrs)	Race	Cause of Death
H0486	Female	49	Caucasian	Anoxia

#### 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 2/27/08



## 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.