

## H2C19.MA / Lot No. 0710456

Human Liver Microsomes

Male, Individual No. 446

0.5 mL at 20 mg protein / mL

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

CYP2C19 Allelic variant	CYP2C19*1/*2
Cytochrome P450 (nmol/mg protein)	0.478
Cytochrome b <sub>5</sub> (nmol/mg protein)	0.539
NADPH-cytochrome c reductase (nmol/mg protein/min)	129 ± 9

Enzyme	Marker substrate reaction (pmol/mg protein/min)	
CYP1A2	Phenacetin <i>O</i> -dealkylation	868 ± 18
CYP2A6	Coumarin 7-hydroxylation	1990 ± 120
CYP2B6	Bupropion hydroxylation	171 ± 8
CYP2C8	Amodiaquine <i>N</i> -dealkylation	2310 ± 30
CYP2C9	Diclofenac 4'-hydroxylation	2180 ± 110
CYP2C19	<i>S</i> -Mephenytoin 4'-hydroxylation	85.5 ± 6.7
CYP2D6	Dextromethorphan <i>O</i> -demethylation	96.3 ± 6.5
CYP2E1	Chlorzoxazone 6-hydroxylation	1120 ± 60
CYP3A4/5	Testosterone 6β-hydroxylation	2780 ± 120
CYP3A4	Midazolam 1'-hydroxylation	293 ± 3
CYP4A11	Lauric acid 12-hydroxylation	2410 ± 70

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

Sample	Gender	Age (yrs)	Race	Cause of Death
H0446	Male	21	Caucasian	Head Trauma

#### 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.