

**H2C9.MA / Lot No. 0710428**

Human Liver Microsomes

Male, Individual No. 539

0.5 mL at 20 mg protein / mL

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

CYP2C9 Allelic variant	CYP2C9*2/*3
Cytochrome P450 (nmol/mg protein)	0.235
Cytochrome b <sub>5</sub> (nmol/mg protein)	0.314
NADPH-cytochrome c reductase (nmol/mg protein/min)	148 ± 6

Enzyme	Marker substrate reaction (pmol/mg protein/min)	
CYP1A2	Phenacetin <i>O</i> -dealkylation	357 ± 6
CYP2A6	Coumarin 7-hydroxylation	314 ± 30
CYP2B6	Bupropion hydroxylation	87.9 ± 4.4
CYP2C8	Amodiaquine <i>N</i> -dealkylation	782 ± 14
CYP2C9	Diclofenac 4'-hydroxylation	1240 ± 170
CYP2C19	<i>S</i> -Mephenytoin 4'-hydroxylation	11.9 ± 0.5
CYP2D6	Dextromethorphan <i>O</i> -demethylation	203 ± 7
CYP2E1	Chlorzoxazone 6-hydroxylation	1440 ± 100
CYP3A4/5	Testosterone 6β-hydroxylation	389 ± 29
CYP3A4	Midazolam 1'-hydroxylation	97.9 ± 9.3
CYP4A11	Lauric acid 12-hydroxylation	1330 ± 50

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

Sample	Gender	Age (yrs)	Race	Cause of Death
H0539	Male	44	Hispanic	Anoxia

**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 2/28/08


**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.