

H2C9.HA / Lot No. 0710439

Human Liver Microsomes

Male, Individual No. 465

0.5 mL at 20 mg protein / mL

Genotype, specific content and activities ^a Content / Rate

CYP2C9 Allelic variant	CYP2C9*1/*1
Cytochrome P450 (nmol/mg protein)	0.436
Cytochrome b ₅ (nmol/mg protein)	0.419
NADPH-cytochrome c reductase (nmol/mg protein/min)	138 ± 1

Enzyme	Marker substrate reaction (pmol/mg protein/min)	
CYP1A2	Phenacetin <i>O</i> -dealkylation	246
CYP2A6	Coumarin 7-hydroxylation	1730 ± 70
CYP2B6	Bupropion hydroxylation	304 ± 17
CYP2C8	Amodiaquine <i>N</i> -dealkylation	2630 ± 90
CYP2C9	Diclofenac 4'-hydroxylation	2960 ± 60
CYP2C19	<i>S</i> -Mephenytoin 4'-hydroxylation	167 ± 8
CYP2D6	Dextromethorphan <i>O</i> -demethylation	166 ± 2
CYP2E1	Chlorzoxazone 6-hydroxylation	730 ± 100
CYP3A4/5	Testosterone 6β-hydroxylation	2570 ± 40
CYP3A4	Midazolam 1'-hydroxylation	516 ± 18
CYP4A11	Lauric acid 12-hydroxylation	1330 ± 30

^a Values for enzyme activities are mean ± standard deviation of three or more determinations.

Sample	Gender	Age (yrs)	Race	Cause of Death
H0465	Male	50	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.