

H2D6.NA / Lot No. 0710445

Human Liver Microsomes

Female, Individual No. 499

0.5 mL at 20 mg protein / mL

Genotype, specific content and activities ^a Content / Rate

CYP2D6 Allelic variant CYP2D6*4/*4

Cytochrome P450 (nmol/mg protein)	0.386
Cytochrome b ₅ (nmol/mg protein)	0.264
NADPH-cytochrome c reductase (nmol/mg protein/min)	118 ± 4

Enzyme	Marker substrate reaction (pmol/mg protein/min)	
CYP1A2	Phenacetin <i>O</i> -dealkylation	308 ± 9
CYP2A6	Coumarin 7-hydroxylation	1180 ± 10
CYP2B6	Bupropion hydroxylation	961 ± 13
CYP2C8	Amodiaquine <i>N</i> -dealkylation	907 ± 87
CYP2C9	Diclofenac 4'-hydroxylation	2530 ± 100
CYP2C19	<i>S</i> -Mephenytoin 4'-hydroxylation	23.0 ± 0.7
CYP2D6	Dextromethorphan <i>O</i> -demethylation	58.9 ± 3.3
CYP2E1	Chlorzoxazone 6-hydroxylation	1260 ± 70
CYP3A4/5	Testosterone 6β-hydroxylation	2120 ± 300
CYP3A4	Midazolam 1'-hydroxylation	107 ± 4
CYP4A11	Lauric acid 12-hydroxylation	1220 ± 40

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

Sample	Gender	Age (yrs)	Race	Cause of Death
H0499	Female	55	Caucasian	Cerebrovascular aneurysm

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 2/27/08



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.