

H2D6.MA / Lot No. 0710425

Human Liver Microsomes
 Female, Individual No. 535
 0.5 mL at 20 mg protein / mL

Genotype, specific content and activities ^a Content / Rate

CYP2D6 Allelic variant	CYP2D6*1/*41
Cytochrome P450 (nmol/mg protein)	0.629
Cytochrome b ₅ (nmol/mg protein)	0.434
NADPH-cytochrome c reductase (nmol/mg protein/min)	150 ± 4

Enzyme	Marker substrate reaction (pmol/mg protein/min)	
CYP1A2	Phenacetin <i>O</i> -dealkylation	1080 ± 70
CYP2A6	Coumarin 7-hydroxylation	804 ± 48
CYP2B6	Bupropion hydroxylation	450 ± 20
CYP2C8	Amodiaquine <i>N</i> -dealkylation	1930 ± 20
CYP2C9	Diclofenac 4'-hydroxylation	2220 ± 80
CYP2C19	<i>S</i> -Mephenytoin 4'-hydroxylation	36.8 ± 1.6
CYP2D6	Dextromethorphan <i>O</i> -demethylation	243 ± 10
CYP2E1	Chlorzoxazone 6-hydroxylation	1740 ± 60
CYP3A4/5	Testosterone 6β-hydroxylation	4280 ± 540
CYP3A4	Midazolam 1'-hydroxylation	367 ± 14
CYP4A11	Lauric acid 12-hydroxylation	1600 ± 30

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

Sample	Gender	Age (yrs)	Race	Cause of Death
H0535	Female	49	Caucasian	Cerebrovascular accident

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.