

H2D6.MA / Lot No. 0710442

Human Liver Microsomes

Male, Individual No. 485

0.5 mL at 20 mg protein / mL

Genotype, specific content and activities ^a Content / Rate

CYP2D6 Allelic variant	CYP2D6*1/*2
Cytochrome P450 (nmol/mg protein)	0.210
Cytochrome b ₅ (nmol/mg protein)	0.224
NADPH-cytochrome c reductase (nmol/mg protein/min)	165 ± 7

Enzyme	Marker substrate reaction (pmol/mg protein/min)	
CYP1A2	Phenacetin <i>O</i> -dealkylation	35.1 ± 4.8
CYP2A6	Coumarin 7-hydroxylation	3.68 ± 0.35
CYP2B6	Bupropion hydroxylation	13.5
CYP2C8	Amodiaquine <i>N</i> -dealkylation	297 ± 22
CYP2C9	Diclofenac 4'-hydroxylation	694
CYP2C19	<i>S</i> -Mephenytoin 4'-hydroxylation	20.4 ± 1.3
CYP2D6	Dextromethorphan <i>O</i> -demethylation	202 ± 4
CYP2E1	Chlorzoxazone 6-hydroxylation	698 ± 164
CYP3A4/5	Testosterone 6β-hydroxylation	42.6 ± 4.0
CYP3A4	Midazolam 1'-hydroxylation	5.07 ± 0.37
CYP4A11	Lauric acid 12-hydroxylation	438 ± 1

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

Sample	Gender	Age (yrs)	Race	Cause of Death
H0485	Male	10	Caucasian	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 2/27/08


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.