

H3A5.NA / Lot No. 0710253

Human Liver Microsomes

Male, Individual No. 307

0.5 mL at 20 mg protein / mL

Genotype, specific content and activities ^a Content / Rate

CYP3A5 Allelic variant	CYP3A5*3/*3
Cytochrome P450 (nmol/mg protein)	0.588
Cytochrome b ₅ (nmol/mg protein)	0.414
NADPH-cytochrome c reductase (nmol/mg protein/min)	316 ± 4

Enzyme	Marker substrate reaction (pmol/mg protein/min)	
CYP1A2	Phenacetin <i>O</i> -dealkylation	606 ± 12
CYP2A6	Coumarin 7-hydroxylation	2160 ± 190
CYP2B6	Bupropion hydroxylation	3400 ± 40
CYP2C8	Amodiaquine <i>N</i> -dealkylation	5150 ± 110
CYP2C9	Diclofenac 4'-hydroxylation	2670 ± 130
CYP2C19	<i>S</i> -Mephenytoin 4'-hydroxylation	29.8 ± 1.5
CYP2D6	Dextromethorphan <i>O</i> -demethylation	279 ± 3
CYP2E1	Chlorzoxazone 6-hydroxylation	12800 ± 900
CYP3A4/5	Testosterone 6β-hydroxylation	4550 ± 440
CYP3A4	Midazolam 1'-hydroxylation	977 ± 41
CYP4A11	Lauric acid 12-hydroxylation	3500 ± 200

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

Sample	Gender	Age (yrs)	Race	Cause of Death
H0307	Male	39	Caucasian	Anoxia

Serology information

- This donor tested negative 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 6/5/07



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.