

## H3A5.HA / Lot No. 0710461

Human Liver Microsomes  
 Female, Individual No. 324  
 0.5 mL at 20 mg protein / mL

### Genotype, specific content and activities <sup>a</sup> Content / Rate

CYP3A5 Allelic variant	CYP3A5*1/*1
Cytochrome P450 (nmol/mg protein)	0.420
Cytochrome b <sub>5</sub> (nmol/mg protein)	0.394
NADPH-cytochrome c reductase (nmol/mg protein/min)	174 ± 1

Enzyme	Marker substrate reaction (pmol/mg protein/min)	
CYP1A2	Phenacetin O-dealkylation	435 ± 11
CYP2A6	Coumarin 7-hydroxylation	270 ± 29
CYP2B6	Bupropion hydroxylation	98.8 ± 3.3
CYP2C8	Amodiaquine N-dealkylation	2160 ± 70
CYP2C9	Diclofenac 4'-hydroxylation	2680 ± 90
CYP2C19	S-Mephenytoin 4'-hydroxylation	122 ± 4
CYP2D6	Dextromethorphan O-demethylation	174 ± 1
CYP2E1	Chlorzoxazone 6-hydroxylation	4670 ± 80
CYP3A4/5	Testosterone 6β-hydroxylation	2050 ± 100
CYP3A4	Midazolam 1'-hydroxylation	383 ± 15
CYP4A11	Lauric acid 12-hydroxylation	889 ± 21

<sup>a</sup> Values for enzyme activities are mean ± standard deviation of three or more determinations.

Sample	Gender	Age (yrs)	Race	Cause of Death
H0324	Female	53	African-American	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 11 June 2013



## 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.