

## HU1A1.NA Lot No. 0710441

Human Liver Microsomes

Male, Individual No. 483

0.5 mL at 20 mg protein / mL

Genotype, specific content and activities <sup>a</sup>		Content / Rate
UGT1A1 Allelic variant		UGT1A1*28/*28
Cytochrome P450 (nmol/mg protein)		0.210
Cytochrome b <sub>5</sub> (nmol/mg protein)		0.269
NADPH-cytochrome c reductase (nmol/mg protein/min)		106 ± 2
Enzyme	Marker substrate reaction (pmol/mg protein/min)	
CYP1A2	Phenacetin <i>O</i> -dealkylation	376 ± 31
CYP2A6	Coumarin 7-hydroxylation	419 ± 7
CYP2B6	Bupropion hydroxylation	32.4 ± 6.3
CYP2C8	Amodiaquine <i>N</i> -dealkylation	383 ± 10
CYP2C9	Diclofenac 4'-hydroxylation	365 ± 26
CYP2C19	<i>S</i> -Mephenytoin 4'-hydroxylation	16.1 ± 0.8
CYP2D6	Dextromethorphan <i>O</i> -demethylation	258 ± 3
CYP2E1	Chlorzoxazone 6-hydroxylation	1140 ± 240
CYP3A4/5	Testosterone 6β-hydroxylation	441 ± 45
CYP3A4	Midazolam 1'-hydroxylation	113 ± 3
CYP4A11	Lauric acid 12-hydroxylation	1080 ± 50
UGT1A1	17β-Estradiol 3-glucuronidation	205 ± 5

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

Sample	Gender	Age (yrs)	Race	Cause of Death
H0483	Male	61	Hispanic	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.



## Store at -80°C

**CAUTION:** This sample should be considered as a potential biohazard and universal precautions should be followed. Intended for in vitro use only.

These data were generated by and are the property of XENOTECH, LLC. These data are not to be reproduced, published or distributed without the express written consent of XENOTECH, LLC

DATASHEET PREPARED 20-APRIL-10