

**H3A5.HA Lot No. 0710238**

Human Liver Microsomes

Male, Individual No. 238

0.5 mL at 20 mg protein / mL

Suspension medium: 250 mM sucrose



<b>Genotype, Specific Content and Activities <sup>a</sup></b>			<b>Content / Rate</b>
CYP3A5 Allelic variant			CYP3A5*1/*1
Cytochrome P450	(nmol/mg protein)		0.138
Cytochrome b <sub>5</sub>	(nmol/mg protein)		0.302
NADPH-cytochrome c reductase	(nmol/mg protein/min)		145 ± 6
<b>Enzyme</b>	<b>Marker Substrate Reaction</b>	<b>Rate (pmol/mg protein/min)</b>	
CYP1A2	7-Ethoxyresorufin O-dealkylation	18.2 ± 0.9	
CYP2A6	Coumarin 7-hydroxylation	NPD	
CYP2B6	S-Mephenytoin N-demethylation	14.6 ± 4.2	
CYP2C8	Paclitaxel (taxol) 6α-hydroxylation	71.5 ± 7.5	
CYP2C9	Diclofenac 4'-hydroxylation	1050 ± 30	
CYP2C19	S-Mephenytoin 4'-hydroxylation	38.0 ± 0.9	
CYP2D6	Dextromethorphan O-demethylation	278 ± 7	
CYP2E1	Chlorzoxazone 6-hydroxylation	1620 ± 20	
CYP3A4/5	Testosterone 6β-hydroxylation	441 ± 14	
CYP4A11	Lauric acid 12-hydroxylation	731 ± 56	

Characterization is performed when the first lot of a product from a given subcellular fraction (e.g., S9) is prepared. Subsequent lots are subject to a verification test only. Values for enzyme activities were determined at a single substrate concentration and are mean ± standard deviation of three or more determinations.

<b>Sample</b>	<b>Gender</b>	<b>Age (yrs)</b>	<b>Race</b>	<b>Cause of Death</b>
H0238	Male	3 months	African American	Cerebrovascular accident

**Serology information**

- This donor tested positive for cytomegalovirus
- This donor tested negative for HIV, HbsAg, and HCV\*
- This donor tested negative for RPR\*\*

\* Antibody to Human Immunodeficiency 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. These data are not to be reproduced, published or distributed without the express written consent of XenoTech.

Datasheet prepared 06 August 2020