

HU1A9.HA Lot No. 0710410

Human Liver Microsomes

Male, Individual No. 504

0.5 mL at 20 mg protein/mL

Suspension medium: 250 mM sucrose

Genotype, Specific Content and Enzyme Activities		Content / Rate	
UGT1A9 Allelic variant		UGT1A9*1/*1	
Cytochrome P450 content	(nmol/mg protein)	0.730	
Cytochrome b ₅ content	(nmol/mg protein)	0.349	
NADPH-cytochrome c reductase	(nmol/mg protein/min)	182 ± 4	
Enzyme	Marker Substrate Reaction	[S] (µM)	Rate (pmol/mg protein/min)
CYP1A2	Phenacetin O-dealkylation	80	455 ± 46
CYP2A6	Coumarin 7-hydroxylation	50	2350 ± 90
CYP2B6	Bupropion hydroxylation	500	825 ± 17
CYP2C8	Amodiaquine N-dealkylation	20	1840 ± 120
CYP2C9	Diclofenac 4'-hydroxylation	100	412 ± 23
CYP2C19	S-Mephenytoin 4'-hydroxylation	400	88.5 ± 6.5
CYP2D6	Dextromethorphan O-demethylation	80	86.8 ± 1.4
CYP2E1	Chlorzoxazone 6-hydroxylation	500	3500 ± 200
CYP3A4/5	Testosterone 6β-hydroxylation	250	12300 ± 400
CYP3A4/5	Midazolam 1'-hydroxylation	30	1480 ± 30
CYP4A11	Lauric acid 12-hydroxylation	100	1480 ± 60
UGT1A9	Propofol glucuronidation	50	12000 ± 400

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.

To measure cytochrome P450 (CYP) activity, liver microsomes (50 µg/mL) were incubated in triplicate at 37 ± 1°C for 10 minutes in potassium phosphate buffer (50 mM, pH 7.4), containing MgCl₂ (3.0 mM), EDTA (1.0 mM), NADP (1.0 mM), glucose-6-phosphate (5.0 mM), glucose-6-phosphate dehydrogenase (1 Unit/mL) and marker substrate, at the final concentrations indicated. Metabolite formation was determined by validated LC-MS/MS methods with deuterated metabolites as internal standards.

To measure UDP-glucuronosyltransferase (UGT) activity, liver microsomes (10 - 250 µg/mL) were incubated in triplicate at 37 ± 1°C for 5 or 10 minutes in Tris-HCl (100 mM, pH 7.7 at 37°C), CHAPS (0.5 mM), EDTA (1.0 mM), MgCl₂ (10 mM), D-saccharic acid 1,4-lactone (100 µM), uridine diphosphate-glucuronic acid (8.0 mM) and marker substrate at the final concentrations indicated.



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 01 July 2014

Donor Information

Sample	Gender	Age (Yrs)	Race	Cause of Death
504	M	60	Caucasian	Cerebrovascular Accident

Serology information

- Cytomegalovirus: donor tested negative.
- RPR*: donor tested negative
- Donor tested negative for HIV, HTLV, HbsAg, and HCV**

* Rapid Plasma Reagin

** Antibody to Human Immunodeficiency Virus, Antibody to Human T Cell Lymphotropic Virus, Hepatitis B Surface Antigen, Antibody to Hepatitis C Virus, respectively.

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.