

HU1A1.NA Lot No. 1810020

Human Liver Microsomes

Male, Individual No. 1042

0.5 mL at 20 mg protein / mL

Suspension medium: 250 mM sucrose

Genotype, Specific Content and Activities ^a

Content / Rate

UGT1A1 Allelic variant

UGT1A1*28/*28

Cytochrome P450

(nmol/mg protein)

0.277

 Cytochrome b₅

(nmol/mg protein)

0.284

NADPH-cytochrome c reductase

(nmol/mg protein/min)

179 ± 5

Enzyme	Marker Substrate Reaction	[S] (μM)	Rate (pmol/mg protein/min)
CYP1A2	Phenacetin O-dealkylation	80	159 ± 7
CYP2A6	Coumarin 7-hydroxylation	50	407 ± 37
CYP2B6	Bupropion hydroxylation	500	88.4 ± 10.5
CYP2C8	Amodiaquine N-dealkylation	20	1190 ± 30
CYP2C9	Diclofenac 4'-hydroxylation	100	2530 ± 70
CYP2C19	S-Mephenytoin 4'-hydroxylation	400	28.2 ± 3.7
CYP2D6	Dextromethorphan O-demethylation	80	147 ± 12
CYP2E1	Chlorzoxazone 6-hydroxylation	500	2090 ± 120
CYP3A4/5	Testosterone 6β-hydroxylation	30	1910 ± 250
CYP3A4	Midazolam 1'-hydroxylation	250	275 ± 17
CYP4A11	Lauric acid 12-hydroxylation	30	1290 ± 30
UGT1A1	17β-Estradiol 3-glucuronidation	100	232 ± 21

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 ± 2°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 ± 2°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 (10.0 mM) and marker substrate at the final concentrations indicated.

Sample	Gender	Age (yrs)	Race	Cause of Death
H1042	Male	51	Caucasian	Cerebrovascular accident

Serology information

- This donor tested negative 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.

This data sheet serves as a Certificate of Analysis and has been approved by Stephanie Helmstetter, Senior Manager.

Signature and Date: Stephanie Helmstetter 15 February 2022