

H1500 Lot No. 2310343

Human Liver Microsomes

Female, Pool of 10

0.5 mL at 20 mg protein / mL

Suspension medium: 250 mM sucrose

Specific Content and Enzyme Activities		Content / Rate	
Cytochrome P450 content	(nmol/mg protein)	0.327	
Cytochrome b ₅ content	(nmol/mg protein)	0.344	
NADPH-cytochrome c reductase	(nmol/mg protein/min)	190 ± 5	
Glucuronidation of 4-Methylumbelliferone	(nmol/mg protein/min)	117 ± 9	
Enzyme	Marker Substrate Reaction	[S] (µM)	Rate (pmol/mg protein/min)
CYP1A2	Phenacetin O-dealkylation	80	480 ± 53
CYP2A6	Coumarin 7-hydroxylation	50	1060 ± 130
CYP2B6	Bupropion hydroxylation	500	487 ± 31
CYP2C8	Amodiaquine N-dealkylation	20	2000 ± 180
CYP2C9	Diclofenac 4'-hydroxylation	100	2180 ± 150
CYP2C19	S-Mephenytoin 4'-hydroxylation	400	78.4
CYP2D6	Dextromethorphan O-demethylation	80	155 ± 16
CYP2E1	Chlorzoxazone 6-hydroxylation	500	2440 ± 210
CYP3A4/5	Testosterone 6β-hydroxylation	250	3600 ± 140
CYP3A4/5	Midazolam 1'-hydroxylation	30	840 ± 76
CYP4A11	Lauric acid 12-hydroxylation	100	1660 ± 0

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. FMO activity was measured under similar conditions except the protein concentration was 1 mg/mL and the buffer was 49 mM Tricine (pH 8.5).

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 (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.

This data sheet serves as a Certificate of Analysis and has been approved by **Stephanie Helmstetter, Assistant Director.**
Signature and Date: Stephanie Helmstetter 29 January 2024

Donor Information

Sample	Gender	Age (Yrs)	Race	Cause of Death
421	F	53	Caucasian	Anoxia
463	F	58	Caucasian	Anoxia
484	F	55	Caucasian	Cerebrovascular accident
517	F	47	Caucasian	Cerebrovascular accident
548	F	61	Caucasian	Anoxia
585	F	44	Caucasian	Anoxia
594	F	51	Caucasian	Anoxia
780	F	48	African American	Cerebrovascular accident
817	F	31	Caucasian	Anoxia
842	F	60	Caucasian	Anoxia

Serology information

- Cytomegalovirus: 6 donors tested positive.
- RPR*: All donors tested negative
- HIV, HTLV, HbsAg, and HCV**: All donors tested negative.

* 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.