

H1500 Lot No. 1210079

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)		0.402
Cytochrome b ₅ content	(nmol/mg protein)		0.269
NADPH-cytochrome c reductase	(nmol/mg protein/min)		130 ± 7
Enzyme	Marker Substrate Reaction	[S] (μM)	Rate (pmol/mg protein/min)
CYP1A2	Phenacetin O-dealkylation	80	440 ± 23
CYP2A6	Coumarin 7-hydroxylation	50	1360 ± 30
CYP2B6	Bupropion hydroxylation	500	515 ± 12
CYP2C8	Amodiaquine N-dealkylation	20	3210 ± 330
CYP2C9	Diclofenac 4'-hydroxylation	100	2370 ± 40
CYP2C19	S-Mephenytoin 4'-hydroxylation	400	86.0 ± 4.2
CYP2D6	Dextromethorphan O-demethylation	80	195 ± 2
CYP2E1	Chlorzoxazone 6-hydroxylation	500	1520 ± 30
CYP3A4/5	Testosterone 6β-hydroxylation	250	3160 ± 130
CYP3A4/5	Midazolam 1'-hydroxylation	30	786 ± 16
CYP4A11	Lauric acid 12-hydroxylation	100	1200 ± 80
FMO	Benzydamine N-oxygenation	500	1110 ± 190
UGT1A1	17β-Estradiol 3-glucuronidation	100	833 ± 18
UGT1A4	Trifluoperazine glucuronidation	25	818 ± 19
UGT1A6	1-Naphthol glucuronidation	500	10900 ± 600
UGT1A9	Propofol glucuronidation	50	4210 ± 90
UGT2B7	Morphine 3-glucuronidation	1000	3720 ± 50

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. 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 ± 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 15 April 2013

Donor Information

Sample	Gender	Age (Yrs)	Race	Cause of Death
90	F	47	Caucasian	Head trauma
132	F	62	Caucasian	Anoxia
158	F	36	African American	Cerebrovascular accident
208	F	78	Caucasian	Head trauma
225	F	72	Caucasian	Cerebrovascular accident
250	F	50	Caucasian	Cerebrovascular accident
251	F	42	Caucasian	Anoxia
259	F	64	Caucasian	Anoxia
265	F	65	Caucasian	Cerebrovascular accident
272	F	31	Caucasian	Cerebrovascular accident

Serology information

- Cytomegalovirus: 5 donors tested positive.
- RPR*: 7 donors tested negative, 1 donor was non-reactive and 2 donors were not determined
- 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.