

Human Liver Microsomes – Pool of 50

Lot No. 2010065

Human Liver Microsomes

Mixed Gender, Pool of 50

Suspension medium: 250 mM sucrose

H0610 0.5 mL at 20 mg/mL

H0620 1.0 mL at 20 mg/mL

H0630 5.0 mL at 20 mg/mL

H0640 50.0 mL at 20 mg/mL

Specific Content and Enzyme Activities		Content / Rate	
Cytochrome P450 content	(nmol/mg protein)	0.453	
Cytochrome b ₅ content	(nmol/mg protein)	0.369	
NADPH-cytochrome c reductase	(nmol/mg protein/min)	172 ± 3	
Enzyme	Marker Substrate Reaction	[S] (µM)	Rate (pmol/mg protein/min)
CYP1A2	Phenacetin O-dealkylation	80	499 ± 33
CYP2A6	Coumarin 7-hydroxylation	50	1300 ± 20
CYP2B6	Bupropion hydroxylation	500	701 ± 88
CYP2C8	Amodiaquine N-dealkylation	20	2600 ± 580
CYP2C9	Diclofenac 4'-hydroxylation	100	3010 ± 480
CYP2C19	S-Mephenytoin 4'-hydroxylation	400	113 ± 2
CYP2D6	Dextromethorphan O-demethylation	80	273 ± 20
CYP2E1	Chlorzoxazone 6-hydroxylation	500	2490 ± 90
CYP2J2	Ebastine hydroxylation	30	365 ± 15
CYP3A4/5	Testosterone 6β-hydroxylation	250	3040 ± 80
CYP3A4/5	Midazolam 1'-hydroxylation	30	948 ± 104
CYP4A11	Lauric acid 12-hydroxylation	100	2040 ± 50
FMO	Benzylamine N-oxygenation	500	962 ± 22
UGT1A1	17β-Estradiol 3-glucuronidation	100	1250 ± 70
UGT1A3	Chenodeoxycholic acid 24-glucuronidation	300	83.1 ± 3.6
UGT1A4	Trifluoperazine glucuronidation	25	735 ± 23
UGT1A6	1-Naphthol glucuronidation	500	14400 ± 1130
UGT1A9	Propofol glucuronidation	50	4200 ± 90
UGT2B7	Morphine 3-glucuronidation	1000	3690 ± 200
UGT2B17	Testosterone 17-glucuronidation	50	482 ± 17

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 (10.0 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 Sekisui XenoTech. These data are not to be reproduced, published or distributed without the express written consent of Sekisui XenoTech.

Datasheet prepared 17 June 2020

Donor Information

Sample	Gender	Age (Yrs)	Race	Cause of Death
342	F	31	Caucasian	Anoxia
397	M	60	Caucasian	Anoxia
409	F	63	Hispanic	Cerebrovascular Accident
415	M	56	Caucasian	Anoxia
422	M	69	Caucasian	Cerebrovascular Accident
424	F	39	Caucasian	Cerebrovascular Accident
445	F	44	Asian	Cerebrovascular Accident
455	F	37	Caucasian	Cerebrovascular Accident
458	M	48	Caucasian	Head Trauma
467	M	33	Caucasian	Anoxia
471	M	49	Caucasian	Anoxia
485	M	10	Caucasian	Anoxia
486	F	49	Caucasian	Anoxia
487	M	48	Caucasian	Head Trauma
490	F	60	Caucasian	Cerebrovascular Accident
500	F	51	Caucasian	Head Trauma
501	F	58	Caucasian	Anoxia
521	M	56	Caucasian	Anoxia
526	M	34	Caucasian	Head Trauma
528	M	60	Caucasian	Head Trauma
529	M	26	Caucasian	Head Trauma
536	F	34	Caucasian	Anoxia
553	M	74	African American	Cerebrovascular Accident
562	M	65	Caucasian	Cerebrovascular Accident
569	M	60	Caucasian	Anoxia
570	M	49	Caucasian	Cerebrovascular Accident
578	F	56	Hispanic	Cerebrovascular Accident
580	M	77	Caucasian	Cerebrovascular Accident
591	F	11	Caucasian	Head Trauma
592	F	83	Caucasian	Head Trauma
593	M	59	Caucasian	Head Trauma
595	M	37	Caucasian	Head Trauma
675	M	5	Caucasian	Anoxia
689	F	5	Caucasian	Head Trauma
723	F	57	Caucasian	Cerebrovascular Accident
725	F	54	Caucasian	Cerebrovascular Accident
751	M	29	Caucasian	Anoxia
779	F	20	Caucasian	Head Trauma
786	F	27	Caucasian	Head Trauma
802	M	33	Caucasian	Cerebrovascular Accident
811	M	16	Caucasian	Head Trauma
816	M	55	Hispanic	Head Trauma
830	M	65	Caucasian	Cerebrovascular Accident
833	M	48	Hispanic	Anoxia
836	M	39	Caucasian	Head Trauma
962	M	47	Caucasian	Anoxia

These data were generated by and are the property of Sekisui XenoTech. These data are not to be reproduced, published or distributed without the express written consent of Sekisui XenoTech.

Donor Information

Sample	Gender	Age (Yrs)	Race	Cause of Death
973	F	63	Caucasian	Cerebrovascular Accident
976	F	61	Caucasian	Head Trauma
997	M	63	Caucasian	Anoxia
998	F	63	Caucasian	Cerebrovascular Accident

Serology information

- Antibody to Cytomegalovirus: 29 of 50 donors tested positive.
- RPR*: 50 donors tested negative.
- HIV, HbsAg, and HCV**: All donors tested negative.

* Rapid Plasma Reagin

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

These data were generated by and are the property of Sekisui XenoTech. These data are not to be reproduced, published or distributed without the express written consent of Sekisui XenoTech.