

H0610.I Lot No. 2210287

Human Intestine Microsomes
Mixed Gender, Pool of 13
150 μ L at 10 mg protein / mL
Suspension medium: 250 mM sucrose

Enzyme Activities		Rate
NADPH-cytochrome c reductase	(nmol/mg protein/min)	31.5 \pm 0.9
Testosterone 6 β -hydroxylation	(pmol/mg protein/min)	810 \pm 21
Midazolam 1'-hydroxylation	(pmol/mg protein/min)	208 \pm 3
Glucuronidation of 4-Methylumbelliferone	(nmol/mg protein/min)	4.36 \pm 0.40

Values for enzyme activities were determined at a single substrate concentration and are mean \pm standard deviation of three or more determinations.

Aprotinin, Leupeptin, and Phenylmethylsulfonyl-fluoride were used in the preparation of these microsomes. Subcellular fractions were prepared from duodenal and jejunal tissue.

To measure cytochrome P450 (CYP) activity, intestine microsomes (0.1 mg/mL) were incubated in triplicate at 37 \pm 2 $^{\circ}$ 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 testosterone (250 μ M), at the final concentrations indicated. Metabolite formation was determined by validated LC-MS/MS methods with deuterated metabolites as internal standards.

To measure cytochrome P450 (CYP) activity, intestine microsomes (0.1 mg/mL) were incubated in triplicate at 37 \pm 2 $^{\circ}$ 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 midazolam (30 μ M), 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, intestine microsomes (0.2 mg/mL) were incubated in triplicate at 37 \pm 2 $^{\circ}$ C for 10 minutes in Tris-HCl (100 mM, pH 7.7 at 37 $^{\circ}$ C), CHAPS (0.5 mM), EDTA (1.0 mM), MgCl₂ (10 mM), D-saccharic acid 1,4-lactone (100 μ M), uridine diphosphate-glucuronic acid (8.0 mM) and 4-methylumbelliferone (1 mM), at the final concentrations indicated. Metabolite formation was determined by validated LC-MS/MS methods with deuterated metabolites as internal standards.



Store at -80 $^{\circ}$ 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 18 November 2022

Donor Information

Sample	Gender	Age (Yrs)	Race	Cause of Death
157	M	57	Caucasian	Anoxia
158	M	58	Caucasian	Head trauma
159	F	49	Caucasian	Cerebrovascular accident
160	M	54	Hispanic	Head trauma
162	M	65	Caucasian	Anoxia
163	M	28	African American	Anoxia
164	F	43	Caucasian	Anoxia
166	F	58	Caucasian	Cerebrovascular accident
167	M	64	Caucasian	Anoxia
168	F	53	African American	Head trauma
169	M	52	Caucasian	Head trauma
170	F	62	Caucasian	Head trauma
171	F	19	Caucasian	Head trauma

Serology information

- Cytomegalovirus: 8 donors tested positive.
- RPR*: All 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 XenoTech. These data are not to be reproduced, published or distributed without the express written consent of XenoTech.