

## H0610.I Lot No. 2310319

Human Intestine Microsomes  
Mixed Gender, Pool of 15  
150  $\mu$ L at 10 mg protein / mL  
Suspension medium: 250 mM sucrose

<b>Enzyme Activities</b>		<b>Rate</b>
NADPH-cytochrome c reductase	(nmol/mg protein/min)	44.5 $\pm$ 3.4
Testosterone 6 $\beta$ -hydroxylation	(pmol/mg protein/min)	1330 $\pm$ 90
Midazolam 1'-hydroxylation	(pmol/mg protein/min)	233 $\pm$ 10

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°C for 10 minutes in potassium phosphate buffer (50 mM, pH 7.4), containing MgCl<sub>2</sub> (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  $\mu$ 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°C for 10 minutes in potassium phosphate buffer (50 mM, pH 7.4), containing MgCl<sub>2</sub> (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  $\mu$ M), at the final concentrations indicated. Metabolite formation was determined by validated LC-MS/MS methods with deuterated metabolites as internal standards.



### 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 22 November 2023

## Donor Information

Sample	Gender	Age (Yrs)	Race	Cause of Death
110	M	70	Caucasian	Cerebrovascular accident
112	M	35	Caucasian	Cerebrovascular accident
113	M	45	Caucasian	Head trauma
114	M	18	Caucasian	Head trauma
115	M	36	Caucasian	Head trauma
172	F	66	Caucasian	Cerebrovascular accident
173	M	57	Caucasian	Cerebrovascular accident
174	F	14	Caucasian	Anoxia
175	M	66	Caucasian	Cerebrovascular accident
176	M	16	Caucasian	Head trauma
177	M	65	Caucasian	Head trauma
178	M	55	Caucasian	Head trauma
179	F	56	Caucasian	Anoxia
180	M	52	Caucasian	Cerebrovascular accident
181	M	55	Caucasian	Cerebrovascular accident

### **Serology information**

- Cytomegalovirus: 9 donors tested positive, 5 donors tested negative, and 1 donor was not determined
- 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.