

H0610.IS9 Lot No. 2210203

Human Intestine S9 Fraction

Mixed Gender, Pool of 39

1.0 mL at 4 mg protein / mL

Suspension medium: 50 mM Tris-HCl, 150 mM KCl, 1 mM EDTA, 20% glycerol, heparin, PMSF, leupeptin, DTT, aprotinin

Enzyme Activities		Rate
NADPH-cytochrome c reductase	(nmol/mg protein/min)	18.8 ± 1.0
Testosterone 6β-hydroxylation	(pmol/mg protein/min)	474 ± 27
Midazolam 1'-hydroxylation	(pmol/mg protein/min)	89.9 ± 8.0
Glucuronidation of 4-Methylumbelliferone	(nmol/mg protein/min)	1.61 ± 0.02

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

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

To measure cytochrome P450 (CYP) activity, intestine S9 samples (0.2 mg/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 testosterone (250 μM), 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 S9 samples (0.2 mg/mL) were incubated in triplicate at 37 ± 2°C for 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 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.

Donor Information

Serology information

- Cytomegalovirus: 25 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.

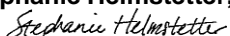


Store at -80°C

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, Senior Manager.

Signature and Date:  26 August 2022

Sample	Gender	Age (Yrs)	Race	Cause of Death
91	F	55	Caucasian	Cerebrovascular accident
94	M	46	Caucasian	Cerebrovascular accident
96	M	51	Caucasian	Head trauma
98	M	27	Caucasian	Head trauma
99	F	44	Caucasian	Cerebrovascular accident
100	F	37	Caucasian	Head trauma
102	M	18	Caucasian	Head trauma
104	F	53	Caucasian	Cerebrovascular accident
105	M	32	Hispanic	Anoxia
106	M	27	Caucasian	Head trauma
107	M	49	Caucasian	Head trauma
135	F	57	Caucasian	Cerebrovascular accident
136	F	57	Caucasian	Head trauma
138	M	54	Caucasian	Head trauma
139	M	63	Hispanic	Cerebrovascular accident
140	F	54	Caucasian	Cerebrovascular accident
142	F	69	Caucasian	Cerebrovascular accident
143	M	42	Caucasian	Cerebrovascular accident
144	M	26	Caucasian	Head trauma
145	M	55	Caucasian	Anoxia
146	M	40	Asian	Anoxia
147	F	47	Caucasian	Head trauma
148	F	61	Caucasian	Cerebrovascular accident
149	F	53	Caucasian	Cerebrovascular accident
151	M	61	Caucasian	Head trauma
152	F	66	Caucasian	Cerebrovascular accident
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

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.