

H0610.ES9 Lot No. 2110214

Human Epidermal Skin S9 Fraction

Mixed Gender, Pool of 6

0.5 mL at 4 mg protein / mL

Suspension medium: 50 mM Tris-HCl, 150 mM KCl, 2 mM EDTA

Enzyme Activities		Rate
NADPH-cytochrome c reductase	(nmol/mg protein/min)	1.10 ± 0.00
Testosterone 6β-hydroxylation	(pmol/mg protein/min)	8.22 ± 0.49
Glucuronidation of 4-Methylumbelliferone	(nmol/mg protein/min)	< 0.1
6α-Methylprednisolone 21-hemisuccinate hydrolysis	(pmol/mg protein/min)	525 ± 49

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, skin 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, skin 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.

To measure carboxylesterase activity, skin S9 samples (0.15 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), and 6α-methylprednisolone 21-hemisuccinate (750 μM), at the final concentrations indicated. Metabolite formation was determined by 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.

Datasheet prepared 11 August 2021

Donor Information

Sample	Gender	Age (Yrs)	Race	Cause of Death
30	F	46	Caucasian	Cerebrovascular accident
32	F	55	Caucasian	Anoxia
33	F	55	Caucasian	Anoxia
34	M	63	Caucasian	Anoxia
35	F	63	Caucasian	Anoxia
36	M	56	Caucasian	Anoxia

Serology information

- Donors were not tested for cytomegalovirus
- All donors tested negative for RPR*
- All donors tested negative for HIV, HbsAg, and HCV**

* 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.