

## H0610.ES9 Lot No. 2310320

Human Epidermal Skin S9 Fraction

Mixed Gender, Pool of 7

0.5 mL at 4 mg protein / mL

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

Enzyme Activities	Rate
6 $\alpha$ -Methylprednisolone 21-hemisuccinate hydrolysis (pmol/mg protein/min)	455 $\pm$ 48

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

To measure carboxylesterase activity, skin S9 samples (0.15 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), and 6 $\alpha$ -methylprednisolone 21-hemisuccinate (750  $\mu$ M), at the final concentrations indicated. Metabolite formation was determined by LC-MS/MS methods with deuterated metabolites as internal standards.

### Donor Information

Sample	Gender	Age (Yrs)	Race	Cause of Death
46	M	35	Caucasian	Anoxia
47	M	68	Caucasian	Anoxia
48	M	66	Caucasian	Anoxia
49	M	46	Caucasian	Cerebrovascular accident
51	F	61	Caucasian	Anoxia
52	F	56	Caucasian	Anoxia
53	F	26	Caucasian	Head trauma

### Serology information

- 1 donor tested negative for cytomegalovirus. 6 donors were not tested.
- 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.



### 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 08 December 2023