

H0610.IS9(NP) Lot No. 1710039

Human Intestine S9 Fraction – PMSF-free

Mixed Gender, Pool of 4

1.0 mL at 4 mg protein / mL

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

| Enzyme Activities | Rate |
|---|-------------|
| 6α-Methylprednisolone 21-hemisuccinate hydrolysis (pmol/mg protein/min) | 4564 ± 114 |

Characterization is performed when the first lot of a product from a given subcellular fraction (e.g., S9) is prepared. Subsequent lots are subject to a verification test only. Values for enzyme activities were determined at a single substrate concentration and are mean ± standard deviation of three or more determinations.

Aprotinin and Leupeptin were used in the preparation of this S9 fraction. Phenylmethylsulfonyl-fluoride was not used in the preparation of this S9 fraction. Subcellular fractions were prepared from duodenal and jejunal tissue.

To measure carboxylesterase activity, intestine S9 samples (0.15 mg/mL) were incubated in triplicate at 37 ± 1°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.

Donor Information

| Sample | Gender | Age (Yrs) | Race | Cause of Death |
|--------|--------|-----------|------------------|--------------------------|
| 153 | M | 21 | African American | Anoxia |
| 154 | F | 49 | Caucasian | Cerebrovascular accident |
| 155 | F | 53 | African American | Cerebrovascular accident |
| 156 | M | 67 | Caucasian | Cerebrovascular accident |

Serology information

- Cytomegalovirus: All 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.

Datasheet prepared 22 May 2017