

R1000.IS9(NP) Lot No. 2310170

Sprague Dawley (SD) Rat Intestine S9 Fraction – PMSF-free
Untreated, Male, Pool of 100

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
6 α -Methylprednisolone 21-hemisuccinate hydrolysis (pmol/mg protein/min)	57870 \pm 1450

Values for enzyme activities were determined at a single substrate concentration and are mean \pm 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 \pm 2°C for 10 minutes in potassium phosphate buffer (50 mM, pH 7.4), containing MgCl₂ (3.0 mM), EDTA (1.0 mM), and 6 α -methylprednisolone 21-hemisuccinate (750 μ M), at the final concentrations indicated. Metabolite formation was determined by LC-MS/MS methods with deuterated metabolites as internal standards.

Animal Information

Species: Rat
Strain: International Genetic Standard (IGS), Sprague Dawley
Sex: Male
Age: ~8 weeks
Vendor: Charles River, Raleigh, NC

Animals were housed in an AAALAC-accredited facility and allowed to acclimate \geq seven days before use.

Food: Purina 5L79 (*ad libitum*)
Water: Automatic watering system (*ad libitum*)
Light/dark cycle: 5:00 am - 5:00 pm, light; 5:00 pm - 5:00 am, dark (12-hour light/dark)
Temperature: 70°F \pm 2°F
Humidity: 30-70 %
Bedding: Beta Chip (hardwood), NEPCO, Warrensburg, NY
Cage: Polycarbonate Shoebox Cage, conventional cage



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 28 June 2023