

R1000.I(NP) Lot No. 2110135

Sprague Dawley (SD) Rat Intestine Microsomes – PMSF-free
Untreated, Male, Pool of 200
150 µL at 10 mg protein / mL
Suspension medium: 250 mM sucrose

Enzyme Activities		Rate
6α-Methylprednisolone 21-hemisuccinate hydrolysis	(pmol/mg protein/min)	20406 ± 549

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 microsomes (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.

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 ≥ 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 ± 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.

Datasheet prepared 18 May 2021