

M1000.I(NP) Lot No. 2310339

CD1 Mouse 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)	98500 ± 4070

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 microsomal fraction. Phenylmethylsulfonyl-fluoride was not used in the preparation of this microsomal 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: Mouse
 Strain: CD1
 Sex: Male
 Age: ~11 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.

This data sheet serves as a Certificate of Analysis and has been approved by Stephanie Helmstetter, Assistant Director.
Signature and Date: Stephanie Helmstetter 15 December 2023