

P2000.I(NP) Lot No. 1510021

Cynomolgus Monkey Intestine Microsomes – PMSF-free
Untreated, Male, Pool of 6
150 µL at 10 mg protein / mL
Suspension medium: 250 mM sucrose

Enzyme Activities	Rate
6α-Methylprednisolone 21-hemisuccinate hydrolysis (pmol/mg protein/min)	4644 ± 755

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 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 ± 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.

Animal Information

Species: Monkey
Strain: Cynomolgus
Sex: Male
Age: 2-3 years
Vendor: Charles River, Reno, NV

Animals were housed in an AAALAC-accredited facility.

Imported animals were quarantined for one month prior to shipment into the United States to reduce the risk of importing Ebola virus-infected monkeys. All animals were under veterinary care and were asymptomatic at the time of euthanasia. All of the monkeys tested negative for Simian Retrovirus. None of the animals examined tested positive for any other infectious agents.



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 14 October 2015