

D1000.I(NP) Lot No. 1310234

Beagle Dog Intestine Microsomes – PMSF-free

Untreated, Male, Pool of 3

150 μ L at 10 mg protein / mL

Suspension medium: 250 mM sucrose

Enzyme Activities	Rate
6 α -Methylprednisolone 21-hemisuccinate hydrolysis (pmol/mg protein/min)	1140 \pm 110

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 \pm standard deviation of three or more determinations.

Aprotinin and Leupeptin were used in the preparation of these microsomes. Phenylmethylsulfonyl-fluoride was not used in the preparation of these microsomes. 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 \pm 1°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: Dog; *Canis familiaris*
 Strain: Beagle
 Sex: Male
 Age: >6 months
 Vendor: Xenometrics, Stilwell, KS

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

Food: Nutrena (*ad libitum*)
 Water: Automatic watering system, tap water (*ad libitum*)
 Light/dark cycle: Not monitored
 Temperature: Ranges from 62°-82°F
 Humidity: Not monitored
 Cage: Indoor/outdoor run cages, plastic coated rod bottom, sanitized at least every 2 weeks



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 26 November 2013