

## D1000.IS9(NP) Lot No. 1310232

Beagle Dog Intestine S9 Fraction – PMSF-free  
Untreated, Male, Pool of 3

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

<i>Enzyme Activities</i>	<i>Rate</i>
6 $\alpha$ -Methylprednisolone 21-hemisuccinate hydrolysis (pmol/mg protein/min)	904 $\pm$ 41

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 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  $\pm$  1°C for 10 minutes in potassium phosphate buffer (50 mM, pH 7.4), containing MgCl<sub>2</sub> (3.0 mM), EDTA (1.0 mM), and 6 $\alpha$ -methylprednisolone 21-hemisuccinate (750  $\mu$ 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