

H0610.I(NP) Lot No. 1410066

Human Intestine Microsomes – PMSF-free
 Mixed Gender, 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)	3860 ± 340

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 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 ± 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.

Donor Information

Sample	Gender	Age (Yrs)	Race	Cause of Death
117	F	53	Caucasian	Cerebrovascular accident
118	F	34	African American	Cerebrovascular accident
119	M	15	Caucasian	Head Trauma
124	M	36	Caucasian	Head Trauma
125	F	66	Caucasian	Cerebrovascular accident
126	M	28	African American	Head Trauma

Serology information

- Cytomegalovirus: 3 donors tested positive.
- RPR*: All donors tested negative.
- HIV, HTLV, HbsAg, and HCV**: All donors tested negative.

* Rapid Plasma Reagin

** Antibody to Human Immunodeficiency Virus, Antibody to Human T Cell Lymphotropic Virus, Hepatitis B Surface Antigen, Antibody to Hepatitis C Virus, respectively.



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 30 May 2014