

**P2000.I Lot No. 1110390**

Cynomolgus Monkey Intestinal Microsomes

Untreated, Male, Pool of 7

150  $\mu$ L at 10 mg protein / mL<sup>b</sup>

Suspension medium: 250 mM sucrose

Specific activities <sup>a</sup>	Rate
NADPH-cytochrome <i>c</i> reductase (nmol/mg protein/min)	81.6 $\pm$ 1.6
Testosterone 6 $\beta$ -hydroxylation (pmol/mg protein/min)	1790 $\pm$ 60
4-Methylumbelliferone glucuronidation (nmol/mg protein/min)	61.3 $\pm$ 1.7

<sup>a</sup> 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.

<sup>b</sup> Aprotinin, Leupeptin, and Phenylmethylsulfonyl-fluoride were used in the preparation of these microsomes. Subcellular fractions were prepared from duodenal and jejunal tissue.

**Animal Information**

Species: Monkey  
Strain: Cynomolgus  
Sex: Male  
Age: 2-6 years  
Vendor: Charles River, Sparks, NV; Charles River, Houston, TX

Monkeys 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, LLC. These data are not to be reproduced, published or distributed without the express written consent of XENOTECH, LLC.

Datasheet prepared 9 January 2012