



**P2000.IS9 / Lot No. 0510273**  
Cynomolgus Monkey Intestinal S9  
Untreated, Male, Pool of 11  
1.0 mL at 4 mg protein / mL<sup>a</sup>

Specific activities <sup>b</sup>	Rate
NADPH-Cytochrome <i>c</i> Reductase (nmol/mg protein/min)	20.4 ± 0.2
Testosterone 6β-hydroxylation (pmol/mg protein/min)	596 ± 47
Glucuronidation of 4-Methylumbelliferone (nmol/mg protein/min)	22.6 ± 0.1

<sup>a</sup> Samples are diluted in a buffer containing the protease inhibitors: Aprotinin, Leupeptin, and Phenylmethylsulfonyl-fluoride. Subcellular fractions were prepared from duodenal and jejunal tissue.

<sup>b</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 are mean ± standard deviation of three or more determinations.

#### Animal Information

Species: Monkey  
Strain: Cynomolgus  
Sex: Male  
Age: 4-8 years  
Vendor: Sierra Biomedical: Sparks, NV

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:** Although strict measures are taken to ensure that intestines obtained from laboratory animals do not harbor infectious diseases, we recommend that all animal products be handled as potential biohazards and universal precautions be followed.

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 expressed written consent of XENOTECH, LLC.

Datasheet prepared 1/3/06