

**P2000.IS9 / Lot No. 0510141**

Cynomolgus Monkey Intestinal S9

Untreated, Male, Pool of 6

1.0 mL at 4 mg protein / mL^a**Specific activities^b****Rate**

| | |
|--|------------|
| NADPH-Cytochrome <i>c</i> Reductase (nmol/mg protein/min) | 47.9 ± 0.9 |
| Testosterone 6β-hydroxylation (pmol/mg protein/min) | 826 ± 31 |
| Glucuronidation of 4-Methylumbelliferone (nmol/mg protein/min) | 16.5 ± 0.2 |

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.

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

^b Values for enzyme activities are mean ± standard deviation of three or more determinations.

**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 7/29/05