

P2000.IS9 Lot No. 1110388

Cynomolgus Monkey Intestinal S9

Untreated, Male, Pool of 7

1.0 mL at 4 mg protein / mL^b

Suspension medium: 50 mM Tris·HCl, 150 mM KCl, 1 mM EDTA, 20% glycerol, heparin, PMSF, leupeptin, DTT, aprotinin

Specific activities^a**Rate**

| | |
|---|------------|
| NADPH-cytochrome <i>c</i> reductase (nmol/mg protein/min) | 46.8 ± 0.2 |
| Testosterone 6β-hydroxylation (pmol/mg protein/min) | 1130 ± 50 |
| 4-Methylumbelliferone glucuronidation (nmol/mg protein/min) | 27.0 ± 0.1 |

^a 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.

^b 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