

P2000.IC Lot No. 1410046

Cynomolgus Monkey Intestine Cytosol

Untreated, Male, Pool of 3

1.0 mL at 4 mg protein / mL

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

Species: Monkey
Strain: Cynomolgus
Sex: Male
Age: 6- 7 years
Vendor: BioCHEMED, Winchester, VA

Animals 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.

Aprotinin, Leupeptin, and Phenylmethylsulfonyl-fluoride were used in the preparation of this cytosol. Subcellular fractions were prepared from duodenal and jejunal tissue.

**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 16 October 2014