

**P2000.C Lot No. 1610269**

Cynomolgus Monkey Liver Cytosol

Untreated, Male, Pool of 4

1.0 mL at 10 mg protein / mL

Suspension medium: 50 mM Tris·HCl, 150 mM KCl, 2 mM EDTA

***Animal Information***

Species: Monkey  
Strain: Cynomolgus  
Sex: Male  
Age: 3-4 years  
Vendor: Worldwide Primates Inc., Miami, FL

Animals were housed in an AAALAC-accredited facility and allowed to acclimate  $\geq$  seven days before use.

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

Datasheet prepared 29 September 2016