

L1000.C Lot No. 1210286

New Zealand Rabbit Liver Cytosol Fraction

Untreated, Male, Pool of 8

1.0 mL at 10 mg protein / mL

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

Animal Information

Species: Rabbit
Strain: New Zealand
Sex: Male
Age: > 9 months
Vendor: Covance Research Products, Denver, PA

Rabbits were laboratory animals and were housed in an AAALAC-accredited facility, which is registered as a research facility with the USDA-APHIS-AC. They were allowed to acclimate for \geq seven days.

Food: Purina 5326 high Fiber Diet (*ad libitum*)
Water: Automatic watering system (*ad libitum*)
Light/dark cycle: 16 hours light / 8 hours dark
Temperature: 68-72 °F
Humidity: 50-55 %
Bedding: None used
Cage: Conventional wire grid cage



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 01 October 2013