

Z6000.C Lot No. 2210264

Gottingen Minipig 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: Miniswine
Strain: Gottingen
Sex: Male
Age: ~ 4 month
Vendor: Marshall BioResources, North Rose, NY

Animals were housed in an AAALAC-accredited facility, which is registered as a research facility with the USDA-APHIS-AC.

Food: Harlan Swine Feed
Water: Automatic watering system (*ad libitum*)
Light/dark cycle: 12-hour light/dark cycle
Temperature: 61-81°F
Humidity: 30-70 %

**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.

This data sheet serves as a Certificate of Analysis and has been approved by **Stephanie Helmstetter, Senior Manager**.
Signature and Date: Stephanie Helmstetter 23 September 2022