

M1000.PC Lot No. 1310042

CD1 Mouse Lung Cytosol

Untreated, Male, Pool of 527

1.0 mL at 5 mg protein / mL

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

Animal Information

Species: Mouse
Strain: CD1
Sex: Male
Age: ~11 weeks
Vendor: Charles River, Raleigh, NC

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

Food: Purina 5L79 (*ad libitum*)
Water: Automatic watering system (*ad libitum*)
Light/dark cycle: 5:00 am - 5:00 pm, light; 5:00 pm - 5:00 am, dark (12-hour light/dark)
Temperature: 70°F \pm 2°F
Humidity: 30-70 %
Cage: Beta Chip (hardwood), NEPCO, Warrensburg, NY

**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 7 August 2013