

R2500.C Lot No. 0410208

Fischer 344 Rat Liver Cytosol

Untreated, Female, Pool of 103

1.0 mL at 10 mg protein / mL

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

Animal Information

Species: Rat
Strain: Fischer 344
Sex: Female
Age: ~ 8 weeks
Vendor: Charles River, Raleigh, NC

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

Food: Harlan Teklad Rodent Chow #8604 (*ad libitum*)
Water: Automatic watering system (*ad libitum*)
Light/dark cycle: 6:00 am - 6:00 pm, light; 6:00 pm - 6:00 am, dark (12-hour light/dark)
Temperature: 72°F \pm 3°F
Humidity: 45-55 %
Bedding: Cell-Sorb Plus (gypsum treated paper product), A&W Products, New Philadelphia, OH
Cage: Polycarbonate Shoebox Cage, conventional 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 02 November 2020