

## R0610.LT Lot No. 2410053

Sprague Dawley (SD) Rat Liver Tritosomes

Tyloxapol-treated, Mixed Gender, Pool of 200

0.25 mL at 2.5 mg protein / mL

Suspension medium: 250 mM Sucrose with 20 mM HEPES, pH 7.4

### Enzyme

Acid Phosphatase	(units <sup>1</sup> /mL)	2.096
Cathepsin B	(units <sup>2</sup> /mg protein)	133 ± 1
RNase Detection		Activity Confirmed

<sup>1</sup> Unit = one unit of acid phosphatase will hydrolyze 1 μmole of 4-nitrophenyl phosphate per minute at pH 4.8 at 37 °C.

<sup>2</sup> Unit = μM AMC released per minute at 38°C.

To measure acid phosphatase content, the following kit was used: Acid phosphatase Assay Kit.  
 Vendor: Sigma-Aldrich. Catalog Number: CS0740.

To detect RNase, the following kit was used: RNaseAlert® Lab Test Kit v2 (modified procedure).  
 Vendor: Life Technologies. Catalog Number: 4479768.

### Animal Information

Species:	Rat	Treatment:	Tyloxapol (Triton WR1339)
Strain:	Sprague Dawley (IGS)	Source:	Sigma (Cat. No. T0307)
Sex:	Mixed Gender	Vehicle:	Sodium Chloride
Age:	~8-9 weeks	Concentration:	100 mg/mL
Vendor:	Charles River, Raleigh, NC	Regimen:	450 mg/kg body weight once on day 1, livers were harvested 120 hours post treatment.

Animals were housed in an AAALAC-accredited facility and allowed to acclimate ≥ 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 ± 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.

This data sheet serves as a Certificate of Analysis and has been approved by **Stephanie Helmstetter, Assistant Director.**  
 Signature and Date: Stephanie Helmstetter 07 March 2024