

## H0610.L Lot No. 1710180

Human Liver Lysosomes

Mixed Gender, Pool of 4

0.25 mL at 2.5 mg protein / mL

Suspension medium: 250 mM Sucrose with 10 mM HEPES, pH 7.0

### Enzyme

Acid Phosphatase	(units/mL)	0.269
Cathepsin B	(units <sup>1</sup> /mg protein)	96.5 ± 3.93
RNase Detection		Activity confirmed

<sup>1</sup> Unit = μM AMC released / min 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.

### Donor Information

Sample	Gender	Age (Yrs)	Race	Cause of Death
1266	F	49	Caucasian	Anoxia
1268	F	53	Caucasian	Anoxia
1271	M	30	African American	Head Trauma
1272	F	20	Caucasian	Head Trauma

#### Serology information

- Antibody to Cytomegalovirus: 3 donors tested positive.
- RPR\*: All donors tested negative.
- HIV(I/II), HbsAg, and HCV\*\*: All donors tested negative.

\* Rapid Plasma Reagin

\*\* Antibody to Human Immunodeficiency Virus, Hepatitis B Surface Antigen, Antibody to Hepatitis C Virus, respectively.



## 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 10 October 2017