

## H1000.C Lot No. 2110107

Human Liver Cytosol

Male, Pool of 10

1.0 mL at 10 mg protein / mL

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

### Donor Information

Sample	Gender	Age (Yrs)	Race	Cause of Death
468	M	47	Caucasian	Cerebrovascular accident
479	M	65	Caucasian	Anoxia
529	M	26	Caucasian	Head Trauma
533	M	28	African American	Anoxia
549	M	55	Caucasian	Cerebrovascular accident
551	M	2	Caucasian	Head Trauma
772	M	46	Caucasian	Anoxia
790	M	52	Caucasian	Cerebrovascular accident
803	M	23	African American	Head Trauma
812	M	73	Caucasian	Cerebrovascular accident

#### Serology information

- Cytomegalovirus: 8 donors tested positive.
- RPR\*: All donors tested negative
- HIV, HTLV, HbsAg, and HCV\*\*: All donors tested negative.

\* Rapid Plasma Reagin

\*\* Antibody to Human Immunodeficiency Virus, Antibody to Human T Cell Lymphotropic 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 17 March 2021