

## H0610.H Lot No. 2310124

Human Liver Homogenate

Mixed Gender, Pool of 20

1.0 mL at 20 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
708	M	56	Caucasian	Head trauma
709	M	50	Caucasian	Cerebrovascular accident
717	M	44	African American	Cerebrovascular accident
718	M	29	Caucasian	Head trauma
720	M	49	Caucasian	Head trauma
721	F	52	Caucasian	Cerebrovascular accident
724	M	61	Caucasian	Cerebrovascular accident
725	F	54	Caucasian	Cerebrovascular accident
727	F	60	Caucasian	Anoxia
732	M	54	Caucasian	Head trauma
733	M	33	Caucasian	Anoxia
734	M	57	Caucasian	Cerebrovascular accident
735	M	34	Caucasian	Anoxia
737	F	49	African American	Anoxia
741	F	74	Caucasian	Head trauma
754	F	66	Caucasian	Cerebrovascular accident
756	F	47	Caucasian	Cerebrovascular accident
759	F	33	Caucasian	Cerebrovascular accident
766	F	38	Caucasian	Anoxia
767	F	24	Caucasian	Cerebrovascular accident

#### Serology information

- Cytomegalovirus: 9 of 20 donors tested positive and 1 donor was not determined
- 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.

This data sheet serves as a Certificate of Analysis and has been approved by **Stephanie Helmstetter, Assistant Director.**

Signature and Date: Stephanie Helmstetter 18 April 2023