

H1500.C / Lot No. 0710245

Human Liver Cytosol

Female, Pool of 10

1.0 mL at 10 mg protein / mL

Human Liver Cytosol

Sample	Gender	Age (Yrs)	Race	Cause of Death
90	F	47	Caucasian	Head Trauma
132	F	62	Caucasian	Anoxia
158	F	36	African American	Cerebrovascular Aneurysm
208	F	78	Caucasian	Head Trauma
225	F	72	Caucasian	Cerebrovascular Aneurysm
250	F	50	Caucasian	Cerebrovascular Stroke
251	F	42	Caucasian	Anoxia
259	F	64	Caucasian	Anoxia
265	F	65	Caucasian	Cerebrovascular Aneurysm
272	F	31	Caucasian	Cerebrovascular Aneurysm

This pool of 10 samples contains liver cytosol from donors 90, 132, 158, 208, 225, 250, 251, 259, 265 and 272. This pool was prepared from identical amounts of liver protein from each donor.

Serology information

- 5 of 10 donors tested positive for cytomegalovirus
- 1 donor tested non-reactive, 2 donors were unknown and 7 donors tested negative for RPR*
- All donors tested negative for HIV, HTLV, HbsAg, and HCV**

* 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***For in vitro use only*

CAUTION: These liver samples are from donors who tested negative for HIV and hepatitis. However, we recommend that these samples be considered as potential biohazards and that universal precautions be used when working with human derived products.

These data were generated by and are the property of XENOTECH, LLC. These data are not to be reproduced, published or distributed without the expressed written consent of XENOTECH, LLC.

DATA SHEET PREPARED 4/6/07