

H1500.C / Lot No. 0610370

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
208	F	78	Caucasian	Head Trauma
231	F	63	Caucasian	Cerebrovascular Aneurysm
239	F	60	Caucasian	Cerebrovascular Aneurysm
256	F	67	Caucasian	Cerebrovascular Aneurysm
324	F	53	African American	Cerebrovascular Aneurysm
338	F	20	Caucasian	Motor Vehicle Accident
340	F	69	Caucasian	Cerebrovascular Aneurysm
353	F	63	Caucasian	Motor Vehicle Accident
356	F	48	Caucasian	Cerebrovascular Aneurysm

This pool of 10 samples contains liver cytosol from donors 90, 208, 231, 239, 256, 324, 338, 340, 353 and 356. This pool was prepared on 6/5/06 from identical amounts of liver protein from each donor.

Serology information

- 3 of 10 donors tested positive for cytomegalovirus and 1 donor was not determined
- 1 donors tested non-reactive, 1 donors were not determined and 8 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 01/17/07