

H1000.C / Lot No. 0710493

Human Liver Cytosol

Male, Pool of 10

1.0 mL at 10 mg protein / mL

Human Liver Cytosol

Sample	Gender	Age (Yrs)	Race	Cause of Death
155	M	46	Hispanic	Head Trauma
314	M	36	Caucasian	Cerebrovascular Aneurysm
315	M	21	Caucasian	Head Trauma
341	M	41	Caucasian	Gun Shot Wound
344	M	58	Caucasian	Anoxia
350	M	33	Hispanic	Head Trauma
351	M	58	Caucasian	Cerebrovascular Aneurysm
357	M	29	Caucasian	Head Trauma
358	M	30	Hispanic	Cerebrovascular Aneurysm
360	M	39	Caucasian	Gun Shot Wound

This pool of 10 samples contains liver cytosol from donors 155, 314, 315, 341, 344, 350, 351, 357, 358 and 360. This pool was prepared from identical amounts of liver protein from each donor.

Serology information

- 6 of 10 donors tested positive for cytomegalovirus
- 0 donors tested non-reactive, 1 donor was not determined and 9 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 12/4/07