

H0610.IC Lot No. 1610315

Human Intestine Cytosol

Mixed Gender, Pool of 15

1.0 mL at 4 mg protein / mL

Suspension medium: 50 mM Tris·HCl, 150 mM KCl, 1 mM EDTA, 20% glycerol, heparin, PMSF, leupeptin, DTT, aprotinin

Donor Information

| Sample | Gender | Age (Yrs) | Race | Cause of Death |
|--------|--------|-----------|-----------|--------------------------|
| 135 | F | 57 | Caucasian | Cerebrovascular accident |
| 136 | F | 57 | Caucasian | Head Trauma |
| 138 | M | 54 | Caucasian | Head Trauma |
| 139 | M | 63 | Hispanic | Cerebrovascular accident |
| 140 | F | 54 | Caucasian | Cerebrovascular accident |
| 142 | F | 69 | Caucasian | Cerebrovascular accident |
| 143 | M | 42 | Caucasian | Cerebrovascular accident |
| 144 | M | 26 | Caucasian | Head Trauma |
| 145 | M | 55 | Caucasian | Anoxia |
| 146 | M | 40 | Asian | Anoxia |
| 147 | F | 47 | Caucasian | Head Trauma |
| 148 | F | 61 | Caucasian | Cerebrovascular accident |
| 149 | F | 53 | Caucasian | Cerebrovascular accident |
| 151 | M | 61 | Caucasian | Head Trauma |
| 152 | F | 66 | Caucasian | Cerebrovascular accident |

Serology information

- Cytomegalovirus: 9 donors tested positive.
- RPR*: All donors tested negative.
- HIV, HbsAg, and HCV**: All donors tested negative.

* Rapid Plasma Reagin

** Antibody to Human Immunodeficiency 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 04 January 2017