

## H0610.IC Lot No. 2010166

Human Intestine Cytosol

Mixed Gender, Pool of 13

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
157	M	57	Caucasian	Anoxia
158	M	58	Caucasian	Head trauma
159	F	49	Caucasian	Cerebrovascular accident
160	M	54	Hispanic	Head trauma
162	M	65	Caucasian	Anoxia
163	M	28	African American	Anoxia
164	F	43	Caucasian	Anoxia
166	F	58	Caucasian	Cerebrovascular accident
167	M	64	Caucasian	Anoxia
168	F	53	African American	Head trauma
169	M	52	Caucasian	Head trauma
170	F	62	Caucasian	Head trauma
171	F	19	Caucasian	Head trauma

#### Serology information

- Cytomegalovirus: 8 of 13 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 13 October 2020