

H0610.IC(NP) Lot No. 1410065

Human Intestine Cytosol – PMSF-free

Mixed Gender, Pool of 6

1.0 mL at 4 mg protein / mL

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

Donor Information

Sample	Gender	Age (Yrs)	Race	Cause of Death
117	F	53	Caucasian	Cerebrovascular accident
118	F	34	African American	Cerebrovascular accident
119	M	15	Caucasian	Head Trauma
124	M	36	Caucasian	Head Trauma
125	F	66	Caucasian	Cerebrovascular accident
126	M	28	African American	Head Trauma

Serology information

- Cytomegalovirus: 3 donors tested positive.
- RPR*: All donors tested negative.
- HIV, HTLV, HbsAg, and HCV**: All donors tested negative.

* 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.

Aprotinin and Leupeptin were used in the preparation of this fraction. Phenylmethylsulfonyl-fluoride was not used in the preparation of this fraction. Subcellular fractions were prepared from duodenal and jejunal tissue.



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 30 May 2014