

H0610.RC Lot No. 1310121

Human Kidney Cytosol

Mixed Gender, Pool of 4

1.0 mL at 5 mg protein / mL

Suspension medium: 50 mM Tris·HCl, 150 mM KCl, 2 mM EDTA

Donor Information

Sample	Gender	Age (Yrs)	Race	Cause of Death
16	M	62	Caucasian	Cerebrovascular accident
18	M	45	Caucasian	Anoxia
21	F	60	Caucasian	Cerebrovascular accident
22	F	60	African American	Anoxia

Serology information

- Cytomegalovirus: Two donors tested positive.
- RPR, 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.

Subcellular fractions were prepared from whole kidney.

**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 5 August 2013