

H0610.PC(S) Lot No. 1410122

Human Lung Cytosol (Smoker)

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	Smoked within past 10 years?
20	M	62	African American	Cerebrovascular Accident	Yes
21	F	27	Caucasian	Cerebrovascular Accident	Yes
27	M	59	Caucasian	Cerebrovascular Accident	Yes
29	F	34	Caucasian	Anoxia	Yes

Serology information

- Cytomegalovirus: 2 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.

**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 23 January 2015