

H0610.PC(NS) Lot No. 1310100

Human Lung Cytosol (Non-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?
9	M	28	Caucasian	Drug overdose	No
14	F	12	Caucasian	Anoxia	No
15	M	65	Caucasian	Cerebrovascular accident	No
17	F	37	Caucasian	Cerebrovascular accident	No

Serology information

- Cytomegalovirus: All donors tested negative
- 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 7 August 2013