

## H0610.PS9(S) Lot No. 1110009

Human Lung S9 (Smoker)

Mixed Gender, Pool of 5

1.0 mL at 5 mg protein / mL

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

Specific activities <sup>a</sup>	Rate
NADPH-cytochrome <i>c</i> reductase (nmol/mg protein/min)	10.6 ± 0.6
7-Ethoxyresorufin <i>O</i> -dealkylation (pmol/mg protein/min)	0.795
Glucuronidation of 4-Methylumbelliferone (nmol/mg protein/min)	< 0.1

<sup>a</sup> Characterization is performed when the first lot of a product from a given subcellular fraction (e.g., S9) is prepared. Subsequent lots are subject to a verification test only. Values for enzyme activities are mean ± standard deviation of three or more determinations.

### Donor Information for Human Lung S9, Lot No. 1110009

Sample	Gender	Age (Yrs)	Race	Cause of Death	Smoked within past 10 years?
2	Female	59	Caucasian	Cerebrovascular accident	Yes
5	Male	19	Caucasian	Head Trauma	Yes
7	Female	18	Caucasian	Head Trauma	Yes
8	Male	40	Caucasian	Cerebrovascular accident	Yes
16	Female	57	Caucasian	Cerebrovascular accident	Yes

#### Serology information

- Two donors tested positive and three donors tested negative for cytomegalovirus
- All donors tested negative for RPR, HIV, HTLV, HbAg, and HCV\*

\* Rapid Plasma Reagin, Antibody to Human Immunodeficiency Virus, Antibody to Human T Cell Lymphotropic Virus, Hepatitis B 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, LLC. These data are not to be reproduced, published or distributed without the express written consent of XENOTECH, LLC.

Datasheet prepared 01 April 2011