

## H0610.P(NS) Lot No. 1310116

Human Lung Microsomes (Non-Smoker)

Mixed Gender, Pool of 4

0.5 mL at 10 mg protein / mL

Suspension medium: 250 mM sucrose

<b>Enzyme Activities</b>		<b>Rate</b>
NADPH-cytochrome c reductase	(nmol/mg protein/min)	27.8 ± 0.8
7-Ethoxyresorufin O-dealkylation	(pmol/mg protein/min)	1.46
Phenacetin O-dealkylation	(pmol/mg protein/min)	<0.5
Glucuronidation of 4-Methylumbelliferone	(nmol/mg protein/min)	<0.1

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 were determined at a single substrate concentration and are mean ± standard deviation of three or more determinations.

To measure cytochrome P450 (CYP) activity, lung microsomes (0.075 mg/mL) were incubated in triplicate at 37 ± 1°C for 10 minutes in potassium phosphate buffer (50 mM, pH 7.4), containing MgCl<sub>2</sub> (3.0 mM), EDTA (1.0 mM), NADP (1.0 mM), glucose-6-phosphate (5.0 mM), glucose-6-phosphate dehydrogenase (1 Unit/mL) and 7-ethoxyresorufin (10 µM), at the final concentrations indicated. Metabolite formation was determined fluorimetrically.

To measure cytochrome P450 (CYP) activity, lung microsomes (0.1 mg/mL) were incubated in triplicate at 37 ± 1°C for 10 minutes in potassium phosphate buffer (50 mM, pH 7.4), containing MgCl<sub>2</sub> (3.0 mM), EDTA (1.0 mM), NADP (1.0 mM), glucose-6-phosphate (5.0 mM), glucose-6-phosphate dehydrogenase (1 Unit/mL) and Phenacetin (80 µM), at the final concentrations indicated. Metabolite formation was determined by validated LC-MS/MS methods with deuterated metabolites as internal standards.

To measure UDP-glucuronosyltransferase (UGT) activity, lung microsomes (0.2 mg/mL) were incubated in triplicate at 37 ± 1°C for 10 minutes in Tris-HCl (100 mM, pH 7.7 at 37°C), CHAPS (0.5 mM), EDTA (1.0 mM), MgCl<sub>2</sub> (10 mM), D-saccharic acid 1,4-lactone (100 µM), uridine diphosphate-glucuronic acid (8.0 mM) and 4-methylumbelliferone (1 mM), at the final concentrations indicated. Metabolite formation was determined by validated LC-MS/MS methods with deuterated metabolites as internal standards.

### Donor Information

<b>Sample</b>	<b>Gender</b>	<b>Age (Yrs)</b>	<b>Race</b>	<b>Cause of Death</b>	<b>Smoked within past 10 years?</b>
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 30 July 2013