

H0610.P(NS) Lot No. 2310312

Human Lung Microsomes (Non-Smoker)

Mixed Gender, Pool of 4

0.5 mL at 10 mg protein / mL

Suspension medium: 250 mM sucrose

| Enzyme Activities | | Rate |
|----------------------------------|-----------------------|-------------|
| NADPH-cytochrome c reductase | (nmol/mg protein/min) | 36.7 ± 0.7 |
| 7-Ethoxyresorufin O-dealkylation | (pmol/mg protein/min) | 0.38 ± 0.08 |
| Phenacetin O-dealkylation | (pmol/mg protein/min) | <2.0 |

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 (0.1 mg/mL) were incubated in triplicate at 37 ± 2°C for 10 minutes in potassium phosphate buffer (50 mM, pH 7.4), containing MgCl₂ (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.

Donor Information

| Sample | Gender | Age (Yrs) | Race | Cause of Death | Smoked within past 10 years? |
|--------|--------|-----------|------------------|----------------|------------------------------|
| 25 | F | 49 | Caucasian | Head Trauma | No |
| 28 | F | 54 | African American | Anoxia | No |
| 39 | M | 65 | Caucasian | Head Trauma | No |
| 41 | M | 57 | Caucasian | Head Trauma | No |

Serology information

- Cytomegalovirus: 2 donors tested positive.
- RPR, HIV, HbsAg, and HCV*: All donors tested negative.

** Rapid Plasma Reagin, Antibody to Human Immunodeficiency 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.

This data sheet serves as a Certificate of Analysis and has been approved by **Stephanie Helmstetter, Assistant Director.**

Signature and Date: Stephanie Helmstetter 07 December 2023