

## M1000 Lot No. 2410002

CD1 Mouse Liver Microsomes Untreated, Male, Pool of 2599 0.5 mL at 20 mg protein / mL

Suspension medium: 250 mM sucrose

| Specific Content and Enzyme Activities  |                       | Content / Rate |
|---|-----------------------|----------------|
| Cytochrome P450 content                 | (nmol/mg protein)     | 0.830          |
| Cytochrome b <sub>5</sub> content       | (nmol/mg protein)     | 0.379          |
| NADPH-cytochrome <i>c</i> reductase     | (nmol/mg protein/min) | 127 ± 4        |
| 7-Ethoxycoumarin <i>O</i> -dealkylation | (pmol/mg protein/min) | 1,970 ± 100    |

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, liver microsomes (50  $\mu$ g/mL) were incubated in triplicate at 37  $\pm$  2°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-ethoxycoumarin (500  $\mu$ M), at the final concentrations indicated. Metabolite formation was determined by validated LC-MS/MS methods with deuterated metabolites as internal standards.

## **Animal Information**

Species: Mouse Strain: CD1 Sex: Male

Age: ~ 11-12 weeks

Vendor: Charles River, Raleigh, NC

Animals were housed in an AAALAC-accredited facility and allowed to acclimate ≥ seven days before use.

Food: Purina 5L79 (ad libitum)

Water: Automatic watering system (ad libitum)

Light/dark cycle: 5:00 am - 5:00 pm, light; 5:00 pm - 5:00 am, dark (12-hour light/dark)

Temperature:  $70^{\circ}\text{F} \pm 2^{\circ}\text{F}$ Humidity:  $30-70^{\circ}$ 

Bedding: Beta Chip (hardwood), NEPCO, Warrensburg, NY Cage: Polycarbonate Shoebox Cage, conventional cage



## 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 29 January 2024