

## R3000 Lot No. 2110132

Wistar Rat Liver Microsomes  
Untreated, Male, Pool of 220  
0.5 mL at 20 mg protein / mL  
Suspension medium: 250 mM sucrose

| <i>Specific Content and Enzyme Activities</i> |                       | <i>Content / Rate</i> |
|---|-----------------------|-----------------------|
| Cytochrome P450 content                       | (nmol/mg protein)     | 0.495                 |
| Cytochrome b <sub>5</sub> content             | (nmol/mg protein)     | 0.464                 |
| NADPH-cytochrome c reductase                  | (nmol/mg protein/min) | 149 ± 10              |
| 7-Ethoxycoumarin O-dealkylation               | (pmol/mg protein/min) | 1620 ± 40             |

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 µg/mL) were incubated in triplicate at 37 ± 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 µM), at the final concentrations indicated. Metabolite formation was determined by validated LC-MS/MS methods with deuterated metabolites as internal standards.

### *Animal Information*

Species: Rat  
Strain: Wistar  
Sex: Male  
Age: ~ 8 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°F ± 2°F  
Humidity: 30-70 %  
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.

Datasheet prepared 23 April 2021