

## M3000 Lot No. 2010026

BALB/c Mouse Liver Microsomes

Untreated, Male, Pool of 398

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)	0.939
Cytochrome b <sub>5</sub> content	(nmol/mg protein)	0.509
NADPH-cytochrome c reductase	(nmol/mg protein/min)	127 ± 8
7-Ethoxycoumarin O-dealkylation	(pmol/mg protein/min)	3770 ± 100

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, 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: Mouse  
 Strain: BALB/c  
 Sex: Male  
 Age: ~ 8-11 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 24 February 2020