

# M1000.S9 / Lot No. 0810364

CD1 Mouse Liver S9

Untreated, Male, Pool of 809

1.0 mL at 20 mg protein / mL

## Specific content and activities<sup>a</sup>

## Content / Rate

Cytochrome P450 (nmol/mg protein)	0.292
Cytochrome b <sub>5</sub> (nmol/mg protein)	0.147
7-Ethoxycoumarin O-dealkylation (pmol/mg protein/min) <sup>b</sup>	680 ± 272
Glucuronidation of 4-methylumbelliferone (nmol/mg protein/min)	77.3 ± 11.2
CDNB (nmol/mg protein/min) <sup>c</sup>	1820 ± 120

<sup>a</sup> 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 are mean ± standard deviation of three or more determinations.

<sup>b</sup> Assay performed at 500 μM substrate concentration.

### Animal Information

Species: Mouse  
 Strain: CD1  
 Sex: Male  
 Age: ~ 11 weeks  
 Vendor: Charles River, Raleigh, NC

Mice were laboratory animals and were housed in an AAALAC-accredited facility, which is registered as a research facility with the USDA-APHIS-AC. They were 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: 45-55 %  
 Bedding: Cell-Sorb Plus (gypsum treated paper product), A&W Products, New Philadelphia, OH  
 Cage: Polycarbonate Shoebox Cage, conventional cage



## Store at -80°C

*For in vitro use only*

**CAUTION:** Although strict measures are taken to ensure that livers obtained from laboratory animals do not harbor infectious diseases, we recommend that all animal products be handled as potential biohazards and universal precautions be followed.

These data were generated by and are the property of XENOTECH, LLC. These data are not to be reproduced, published or distributed without the expressed written consent of XENOTECH, LLC.

Data sheet prepared 28 October 2008