

M1500 Lot No. 1010110

CD1 Mouse Liver Microsomes

Untreated, Female, Pool of 800

0.5 mL at 20 mg protein / mL

Suspension medium: 250 mM sucrose

Specific content and activities ^a	Content / Rate
Cytochrome P450 (nmol/mg protein)	0.872
Cytochrome b ₅ (nmol/mg protein)	0.514
NADPH-cytochrome <i>c</i> reductase (nmol/mg protein/min)	183 ± 3
7-Ethoxycoumarin <i>O</i> -dealkylation (pmol/mg protein/min) ^b	3660 ± 90

^a 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.

^b Assay performed at 500 µM substrate concentration.

Animal Information

Species: Mouse
Strain: CD1
Sex: Female
Age: ~ 9-12 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°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, LLC. These data are not to be reproduced, published or distributed without the express written consent of XENOTECH, LLC.

DATASHEET PREPARED 30-APRIL-10