

R2000 / Lot No. 0610244

Fisher 344 Rat Liver Microsomes

Untreated, Male, Pool of 105

0.5 mL at 20 mg protein / mL

Specific content and activities^a

Content / Rate

Cytochrome P450 (nmol/mg protein)	0.587
Cytochrome b ₅ (nmol/mg protein)	0.479
NADPH-cytochrome <i>c</i> reductase (nmol/mg protein/min)	154 ± 4
7-Ethoxycoumarin <i>O</i> -dealkylation (pmol/mg protein/min) ^b	1190 ± 40

^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 are mean ± standard deviation of three or more determinations.

^b Assay performed at 500 µM substrate concentration. Previous lots were characterized at 100 µM.

Animal Information

Species: Rat
 Strain: Fischer 344
 Sex: Male
 Age: 7-11 weeks
 Vendor: Charles River, Raleigh, NC

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

Food: Harlan Teklad Rodent Chow #8604 (*ad libitum*)
 Water: Automatic watering system (*ad libitum*)
 Light/dark cycle: 6:00 am - 6:00 pm, light; 6:00 pm - 6:00 am, dark (12-hour light/dark)
 Temperature: 72°F ± 3°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 9/8/06