

**R3500 / Lot No. 0010183**

Wistar Rat Liver Microsomes
Untreated, Female, Pool of 50
0.5 mL at 20 mg protein / mL

Specific content and activities ^a	Content / Rate
Cytochrome P450 (nmol/mg protein)	0.436
Cytochrome b ₅ (nmol/mg protein)	0.576
NADPH-cytochrome <i>c</i> reductase (nmol/mg protein/min)	335 ± 3
7-Ethoxycoumarin <i>O</i> -dealkylation (pmol/mg protein/min) ^b	1140 ± 20

^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. Lots prepared prior to 1999 were characterized at 100 µM substrate concentration.

Animal Information

Species: Rat
Strain: Wistar
Sex: Female
Age: ~8 weeks
Vendor: Charles River, Wilmington, MA

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**

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.

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 expressed written consent of XENOTECH, LLC.