

R3000 Lot No. 0910241

Wistar Rat Liver Microsomes

Untreated, Male, Pool of 200

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)	0.780
Cytochrome b ₅	(nmol/mg protein)	0.519
NADPH-cytochrome <i>c</i> reductase	(nmol/mg protein/min)	172 ± 3
7-Ethoxycoumarin <i>O</i> -dealkylation	(pmol/mg protein/min) ^b	2000 ± 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: Rat
 Strain: Wistar
 Sex: Male
 Age: ~ 8 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 16-OCTOBER-09