


R3000.S9 / Lot No. 0510017

Wistar Rat Liver S9

Untreated, Male, Pool of 100

1.0 mL at 20 mg protein / mL

Specific content and activities	Content / Rate
Cytochrome P450 (nmol/mg protein)	0.181
Cytochrome b ₅ (nmol/mg protein)	0.131
7-Ethoxycoumarin <i>O</i> -dealkylation (pmol/mg protein/min) ^a	388 ± 59
Glucuronidation of 4-methylumbelliferone (nmol/mg protein/min)	70.3 ± 3.2
CDNB (nmol/mg protein/min) ^b	623 ± 13

Values for enzyme activities are mean ± standard deviation of three or more determinations.

^a Assay performed at 500 μM substrate concentration.

^b 1-chloro-2,4-dinitrobenzene-glutathione conjugation by Glutathione Transferase

Animal Information

Species: Rat
 Strain: Wistar
 Sex: Male
 Age: ~8 weeks
 Vendor: Charles River, NC

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

Food: Harlan Teklad Rodent Laboratory 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.

DATA SHEET PREPARED 4/5/05