

R3000.S9 / Lot No. 0710487

Wistar Rat Liver S9

Untreated, Male, Pool of 100

1.0 mL at 20 mg protein / mL

Specific content and activities^a

Content / Rate

Cytochrome P450 (nmol/mg protein)	0.172
Cytochrome b ₅ (nmol/mg protein)	0.139
7-Ethoxycoumarin <i>O</i> -dealkylation (pmol/mg protein/min)	400 ± 33
Glucuronidation of 4-methylumbelliferone (nmol/mg protein/min)	77.7 ± 1.2
CDNB (nmol/mg protein/min) ^b	615 ± 10

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

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: 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 10/30/07