

**R1500.S9 Lot No. 1110039**

Sprague Dawley (SD) Rat Liver S9  
Untreated, Female, Pool of 200  
1.0 mL at 20 mg protein / mL  
Suspension medium: 50 mM Tris·HCl,  
150 mM KCl,  
2 mM EDTA

Specific content and activities <sup>a</sup>	Content / Rate
Cytochrome P450 (nmol/mg protein)	0.199
Cytochrome b <sub>5</sub> (nmol/mg protein)	0.150
7-Ethoxycoumarin <i>O</i> -dealkylation (pmol/mg protein/min) <sup>b</sup>	196 ± 6
Glucuronidation of 4-methylumbelliferone (nmol/mg protein/min)	61.6 ± 3.3
CDNB (nmol/mg protein/min) <sup>c</sup>	645 ± 6

<sup>a</sup> Values for enzyme activities are mean ± standard deviation of three or more determinations.

<sup>b</sup> Assay performed at 500 µM substrate concentration.

<sup>c</sup> 1-chloro-2,4-dinitrobenzene-glutathione conjugation by Glutathione Transferase.

**Animal Information**

Species: Rat  
Strain: International Genetic Standard (IGS), Sprague Dawley  
Sex: Female  
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 14 April 11