

**R1000.P / Lot No. 1110008**

Sprague Dawley (SD) Rat Lung Microsomes

Untreated, Male, Pool of 299

0.5 mL at 10 mg protein / mL

Suspension medium: 250 mM sucrose

Specific activities <sup>a</sup>	Rate
NADPH-cytochrome <i>c</i> reductase (nmol/mg protein/min)	34.9 ± 1.3
7-Ethoxyresorufin <i>O</i> -dealkylation (pmol/mg protein/min)	2.17
Glucuronidation of 4-Methylumbelliferone (nmol/mg protein/min)	13.6 ± 0.2

<sup>a</sup> 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.

**Animal Information**

Species: Rat  
Strain: International Genetic Standard (IGS), Sprague Dawley  
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 February 2011