

R1000.RS9 / Lot No. 1010154

Sprague Dawley (SD) Rat Kidney S9

Untreated, Male, Pool of 149

1.0 mL at 5 mg protein / mL

Suspension medium: 50 mM Tris·HCl,
150 mM KCl,
2mM EDTA

Specific activities ^a	Rate
NADPH-cytochrome c reductase (nmol/mg protein/min)	11.4 ± 0.2
Lauric Acid 12-hydroxylation (pmol/mg protein/min)	486 ± 71
Glucuronidation of 4-Methylumbelliferone (nmol/mg protein/min)	26.8 ± 0.6

^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.

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: Harlan Teklad Rodent 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: 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 15 December 2010