

R1000.E / Lot No. 1110137

Sprague Dawley (SD) Rat Skin Microsomes
 Untreated, Male, Pool of 50
 250 µL at 10 mg protein / mL
 Suspension medium: 250 mM sucrose

Specific activities ^a	Rate
NADPH-cytochrome c reductase (nmol/mg protein/min)	17.4 ± 0.0
Testosterone 6β-hydroxylation (pmol/mg protein/min)	9.70 ± 0.59
4-Methylumbelliferone glucuronidation (nmol/mg protein/min)	31.8 ± 0.1
Clopidogrel hydrolysis (pmol/mg protein/min)	193 ± 9
Methylprednisolone 21-hemisuccinate hydrolysis (pmol/mg protein/min)	919 ± 19

^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 were determined at a single substrate concentration and are mean ± standard deviation of three or more determinations.

Subcellular fractions were prepared from dorsal full-thickness skin.

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 30 March 2012