

**M1000.ES9 / Lot No. 1110114**

CD1 Mouse Skin S9 Fraction

Untreated, Male, Pool of 100

1.0 mL at 5 mg protein / mL

Suspension medium: 50 mM Tris-HCl,  
150 mM KCl,  
2 mM EDTA**Specific activities <sup>a</sup>****Rate**

NADPH-cytochrome <i>c</i> reductase (nmol/mg protein/min)	7.11 ± 0.42
Testosterone 6β-hydroxylation (pmol/mg protein/min)	5.83
4-Methylumbelliferone glucuronidation (nmol/mg protein/min)	2.34 ± 0.09
Clopidogrel hydrolysis (pmol/mg protein/min)	46.5 ± 8.6
Methylprednisolone 21-hemisuccinate hydrolysis (pmol/mg protein/min)	842 ± 26

<sup>a</sup> 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: Mouse  
Strain: CD1  
Sex: Male  
Age: ~ 11-12 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