

## M1000.E Lot No. 1810051

CD1 Mouse Skin Microsomes  
 Untreated, Male, Pool of 91  
 250  $\mu$ L at 10 mg protein / mL  
 Suspension medium: 250 mM sucrose

Enzyme Activities		Rate
NADPH-cytochrome c reductase	(nmol/mg protein/min)	11.8 $\pm$ 0.4
Testosterone 6 $\beta$ -hydroxylation	(pmol/mg protein/min)	24.6 $\pm$ 1.9
4-Methylumbelliferone glucuronidation	(nmol/mg protein/min)	15.6 $\pm$ 1.3
Clopidogrel hydrolysis	(pmol/mg protein/min)	41.4 $\pm$ 3.4
Methylprednisolone 21-hemisuccinate hydrolysis	(pmol/mg protein/min)	1006 $\pm$ 15

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  $\pm$  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  $\geq$  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  $\pm$  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. These data are not to be reproduced, published or distributed without the express written consent of XenoTech.

Datasheet prepared 22 May 2018