

Z0610.ES9 Lot No. 1710032

Minipig Skin S9 Fraction

Untreated, Mixed Gender, Pool of 3

0.5 mL at 4 mg protein / mL

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

Specific activities ^a	Rate
NADPH-cytochrome c reductase (nmol/mg protein/min)	2.12 ± 0.12
Testosterone 6β-hydroxylation (pmol/mg protein/min)	4.87 ± 0.31
4-Methylumbelliferone glucuronidation (nmol/mg protein/min)	1.22 ± 0.38
Clopidogrel hydrolysis (pmol/mg protein/min)	20.1 ± 4.5
Methylprednisolone 21-hemisuccinate hydrolysis (pmol/mg protein/min)	625 ± 10

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

Animal Information

Species: Miniswine
 Strain: Gottingen
 Sex: Male (1) Female (2)
 Age: Sexually Mature
 Vendor: Xenometrics, Stilwell, KS

Animals were housed in an AAALAC-accredited facility and allowed to acclimate ≥ seven days before use.

Food: Harlan Swine Feed
 Water: Automatic watering system (*ad libitum*)
 Light/dark cycle: 12 hour light cycle
 Temperature: Ranges from 61°-81°F
 Humidity: 30-70%



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 28 February 2017