

**Z2000 / Lot No. 0310183**

Sinclair Minipig Liver Microsomes

Untreated, Male, Pool of 3

0.5 mL at 20 mg protein / mL

**Specific content and activities<sup>a</sup>****Content / Rate**

Cytochrome P450 (nmol/mg protein)

0.775

Cytochrome b<sub>5</sub> (nmol/mg protein)

0.478

NADPH-cytochrome *c* reductase (nmol/mg protein/min)

161 ± 7

7-Ethoxycoumarin *O*-dealkylation (pmol/mg protein/min)<sup>b</sup>

1190 ± 30

<sup>a</sup> 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.

<sup>b</sup> Assay performed at 500 µM substrate concentration. Lots prepared prior to 1999 were characterized at 100 µM substrate concentration.

**Animal Information**

Species: Miniswine

Breed: Sinclair

Sex: Male

Age: ~3 years old

Vendor: Sinclair research Center, Columbia, MO.

Minipigs were laboratory animals and were housed in an AAALAC-accredited facility, which is registered as a research facility with the USDA-APHIS-AC.

Food: Sinclair S-9 Hog Ration

Water: Tap water (*ad libitum*)

Light/dark cycle: 12 hour light cycle, 6:00 am to 6:00 pm

Temperature: Ranges from 62°-85°F

Humidity: Not monitored

**Store at -80°C***For in vitro use only*

**CAUTION:** Although strict measures are taken to ensure that livers obtained from laboratory animals do not harbor infectious diseases, we recommend that all animal products be handled as potential biohazards and universal precautions be followed.

These data were generated by and are the property of XENOTECH, LLC. These data are not to be reproduced, published or distributed without the expressed written consent of XENOTECH, LLC.

DATA SHEET PREPARED 10/8/03