

**G1000 / Lot No. 0510021**

Hartley Albino Guinea Pig Liver Microsomes
Untreated, Male, Pool of 50
0.5 mL at 20 mg protein / mL

Specific content and activities ^a	Content / Rate
Cytochrome P450 (nmol/mg protein)	0.945
Cytochrome b ₅ (nmol/mg protein)	0.778
NADPH-cytochrome <i>c</i> reductase (nmol/mg protein/min)	185 ± 12
7-Ethoxycoumarin <i>O</i> -dealkylation (pmol/mg protein/min) ^b	1940 ± 60

^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 are mean ± standard deviation of three or more determinations.

^b Assay performed at 500 µM substrate concentration. Lots prepared prior to 1999 were characterized at 100 µM substrate concentration.

Animal Information

Species: Guinea Pig
Strain: Hartley Albino
Sex: Male
Age: ~9 weeks
Vendor: Charles River, NC

Guinea pigs were laboratory animals and were housed in an AAALAC-accredited facility, which is registered as a research facility with the USDA-APHIS-AC. They were allowed to acclimate ≥ seven days before use.

Food: Purina Rodent laboratory Chow #8604 (*ad libitum*)
Water: Automatic watering system (*ad libitum*)
Light/dark cycle: 6:00 am – 6:00 pm light, 6:00 pm – 6:00 am dark (12-hour light/dark)
Temperature: 72°F ± 3°F
Humidity: 45-55%
Bedding: Cell-Sorb Plus (gypsum treated paper product), A&W Products, New Philadelphia, OH
Cage: Polycarbonate Shoebox Cage, conventional cage

**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 3/22/05