

G1000 Lot No. 0910275

Hartley Albino Guinea Pig Liver Microsomes

Untreated, Male, Pool of 37

0.5 mL at 20 mg protein / mL

Suspension medium: 250 mM sucrose

Specific content and activities^a**Content / Rate**

Cytochrome P450	(nmol/mg protein)	0.805
Cytochrome b ₅	(nmol/mg protein)	0.633
NADPH-cytochrome <i>c</i> reductase	(nmol/mg protein/min)	133 ± 3
7-Ethoxycoumarin <i>O</i> -dealkylation	(pmol/mg protein/min) ^b	2410 ± 240

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

^b Assay performed at 500 µM substrate concentration.

Animal Information

Species: Guinea Pig
Strain: Hartley Albino
Sex: Male
Age: ~ 9-10 weeks
Vendor: Charles River, Raleigh, NC

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

Food: Purina 5L97 w/ Vit. C (*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 14-OCTOBER-09