

S1000.S9 / Lot No. 0710394

Golden Syrian Hamster Liver S9

Untreated, Male, Pool of 100

1.0 mL at 20 mg protein / mL

Specific content and activities

Content / Rate

Cytochrome P450 (nmol/mg protein)	0.239
Cytochrome b ₅ (nmol/mg protein)	0.0835
7-Ethoxycoumarin <i>O</i> -dealkylation (pmol/mg protein/min) ^a	1350 ± 30
Glucuronidation of 4-methylumbelliferone (nmol/mg protein/min)	63.0 ± 3.2
CDNB (nmol/mg protein/min) ^b	4270 ± 40

Values for enzyme activities are mean ± standard deviation of three or more determinations.

^a Assay performed at 500 μM substrate concentration.

^b 1-chloro-2,4-dinitrobenzene-glutathione conjugation by Glutathione Transferase

Animal Information

Species: Hamster
 Strain: Golden Syrian
 Sex: Male
 Age: ~8 weeks
 Vendor: Charles River, Raleigh, NC

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

Food: Purina 5L79 (*ad libitum*)
 Water: Automatic watering system (*ad libitum*)
 Light/dark cycle: 6:00 am - 8:00 pm, light; 8:00 pm - 6:00 am, dark (14/10-hour light/dark)
 Temperature: 70°F ± 2°F
 Humidity: 30-70 %
 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 09/27/07