

S1000 Lot No. 2110141

Golden Syrian Hamster Liver Microsomes
Untreated, Male, Pool of 75
0.5 mL at 20 mg protein / mL
Suspension medium: 250 mM sucrose

<i>Specific Content and Enzyme Activities</i>		<i>Content / Rate</i>
Cytochrome P450 content	(nmol/mg protein)	0.822
Cytochrome b ₅ content	(nmol/mg protein)	0.484
NADPH-cytochrome c reductase	(nmol/mg protein/min)	183 ± 6
7-Ethoxycoumarin O-dealkylation	(pmol/mg protein/min)	4990 ± 220

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

To measure cytochrome P450 (CYP) activity, liver microsomes (50 µg/mL) were incubated in triplicate at 37 ± 2°C for 10 minutes in potassium phosphate buffer (50 mM, pH 7.4), containing MgCl₂ (3.0 mM), EDTA (1.0 mM), NADP (1.0 mM), glucose-6-phosphate (5.0 mM), glucose-6-phosphate dehydrogenase (1 Unit/mL) and 7-ethoxycoumarin (500 µM), at the final concentrations indicated. Metabolite formation was determined by validated LC-MS/MS methods with deuterated metabolites as internal standards.

Animal Information

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

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: 14 hours light / 10 hours 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. These data are not to be reproduced, published or distributed without the express written consent of XenoTech.

Datasheet prepared 27 April 2021