

Z6000 Lot No. 2310311

Gottingen Minipig Liver Microsomes Untreated, Male, Pool of 4 0.5 mL at 20 mg protein / mL

Suspension medium: 250 mM sucrose

Specific Content and Enzyme Activities		Content / Rate
Cytochrome P450 content Cytochrome b ₅ content	(nmol/mg protein) (nmol/mg protein)	0.964 0.469
NADPH-cytochrome <i>c</i> reductase 7-Ethoxycoumarin <i>O</i> -dealkylation	(nmol/mg protein/min) (pmol/mg protein/min)	178 ± 4 2670 ± 230

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 μ g/mL) were incubated in triplicate at 37 \pm 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 μ M), at the final concentrations indicated. Metabolite formation was determined by validated LC-MS/MS methods with deuterated metabolites as internal standards.

Animal Information

Species: Miniswine
Strain: Gottingen
Sex: Male
Age: ~ 4 months

Vendor: Marshall BioResources, North Rose, NY

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: Harlan Swine Feed

Water: Automatic watering system (ad libitum)

Light/dark cycle: 12-hour light/dark cycle

Temperature: 61-81°F Humidity: 30-70 %



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.