

Z6000 Lot No. 2210262

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	(nmol/mg protein)	0.746
Cytochrome b ₅ content	(nmol/mg protein)	0.578
NADPH-cytochrome c reductase	(nmol/mg protein/min)	160 ± 4
7-Ethoxycoumarin O-dealkylation	(pmol/mg protein/min)	2560 ± 140

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: 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.

This data sheet serves as a Certificate of Analysis and has been approved by Stephanie Helmstetter, Senior Manager.

Signature and Date: Stephanie Helmstetter 23 September 2022