

L1000 Lot No. 2310309

New Zealand Rabbit Liver Microsomes
 Untreated, Male, Pool of 4
 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)	1.342
Cytochrome b ₅ content	(nmol/mg protein)	0.918
NADPH-cytochrome c reductase	(nmol/mg protein/min)	258 ± 4
7-Ethoxycoumarin O-dealkylation	(pmol/mg protein/min)	2870 ± 240

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: Rabbit
 Strain: New Zealand
 Sex: Male
 Age: >9 months
 Vendor: BioChemed, Winchester, VA

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

Food: Purina 5326 High Fiber Diet (*ad libitum*)
 Water: Automatic watering system (*ad libitum*)
 Light/dark cycle: 16 hours light / 8 hours dark
 Temperature: 68°F -72°F
 Humidity: 50-55 %
 Bedding: None used
 Cage: Conventional wire grid 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.

This data sheet serves as a Certificate of Analysis and has been approved by **Stephanie Helmstetter, Assistant Director.**
 Signature and Date: Stephanie Helmstetter 17 November 2023