

P2000 Lot No. 1810093

Cynomolgus Monkey Liver Microsomes

Untreated, Male, Pool of 6

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)	0.772
Cytochrome b ₅ content	(nmol/mg protein)	0.399
NADPH-cytochrome c reductase	(nmol/mg protein/min)	108 ± 2
7-Ethoxycoumarin O-dealkylation	(pmol/mg protein/min)	3710 ± 50

Characterization is performed when the first lot of a product from a given subcellular fraction (e.g., S9) is prepared. Subsequent lots are subject to a verification test only. 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 ± 1°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: Monkey
 Strain: Cynomolgus
 Sex: Male
 Age: 5-7 years
 Vendor: Xenometrics, Stilwell, KS

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

Imported animals were quarantined for one month prior to shipment into the United States to reduce the risk of importing Ebola virus-infected monkeys. All animals were under veterinary care and were asymptomatic at the time of euthanasia. All of the monkeys tested negative for Simian Retrovirus. None of the animals examined tested positive for any other infectious agents



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 22 May 2018