

P1500.S9 Lot No. 1110079

Rhesus Monkey Liver S9

Untreated, Female, Pool of 5

1.0 mL at 20 mg protein / mL

Suspension medium: 50 mM Tris·HCl,
150 mM KCl,
2 mM EDTA

Specific content and activities ^a	Content / Rate
Cytochrome P450 (nmol/mg protein)	0.333
Cytochrome b ₅ (nmol/mg protein)	0.081
7-Ethoxycoumarin O-dealkylation (pmol/mg protein/min) ^b	993 ± 18
Glucuronidation of 4-methylumbelliferone (nmol/mg protein/min)	81.9 ± 7.3
CDNB (nmol/mg protein/min) ^c	1180 ± 20

^a Values for enzyme activities are mean ± standard deviation of three or more determinations.

^b Assay performed at 500 µM substrate concentration.

^c 1-chloro-2,4-dinitrobenzene-glutathione conjugation by Glutathione Transferase.

Animal Information

Species: Monkey
Strain: Rhesus
Sex: Female
Age: 3-11 years
Vendor: Covance, Alice, TX

Monkeys were housed in an AAALAC-accredited facility

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, LLC. These data are not to be reproduced, published or distributed without the express written consent of XENOTECH, LLC.

Datasheet prepared 01 June 11