

L1500.S9 Lot No. 1010271

New Zealand Rabbit Liver S9

Untreated, Female, Pool of 4

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)	0.533
Cytochrome b ₅ (nmol/mg protein)	0.290
7-Ethoxycoumarin O-dealkylation (pmol/mg protein/min) ^b	693 ± 38
Glucuronidation of 4-methylumbelliferone (nmol/mg protein/min)	66.0 ± 3.9
CDNB (nmol/mg protein/min) ^c	3110 ± 110

^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: Rabbit
Strain: New Zealand
Sex: Female
Age: Sexually Mature
Vendor: Covance Research Products, Denver, PA

Rabbits were laboratory animals and were housed in an AAALAC-accredited facility, which is registered as a research facility with the USDA-APHIS-AC. They were allowed to acclimate for ≥ 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, LLC. These data are not to be reproduced, published or distributed without the express written consent of XENOTECH, LLC.

Datasheet prepared 16 February 11