

L1000.S9 / Lot No. 0710108

New Zealand Rabbit Liver S9

Untreated, Male, Pool of 8

1.0 mL at 20 mg protein / mL

Specific content and activities^a

Content / Rate

Cytochrome P450 (nmol/mg protein)	0.465
Cytochrome b ₅ (nmol/mg protein)	0.339
7-Ethoxycoumarin <i>O</i> -dealkylation (pmol/mg protein/min) ^b	830 ± 21
Glucuronidation of 4-methylumbelliferone (nmol/mg protein/min)	63.5 ± 0.8
CDNB (nmol/mg protein/min) ^c	4450 ± 150

^a 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 are mean ± standard deviation of three or more determinations.

^b Assay performed at 500 µM substrate concentration. Previous lots were characterized at 100 µM.

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

Animal Information

Species: Rabbit
 Strain: New Zealand
 Sex: Male
 Age: > 6 Months
 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

For in vitro use only

CAUTION: Although strict measures are taken to ensure that livers obtained from laboratory animals do not harbor infectious diseases, we recommend that all animal products be handled as potential biohazards and universal precautions be followed.

These data were generated by and are the property of XENOTECH, LLC. These data are not to be reproduced, published or distributed without the expressed written consent of XENOTECH, LLC.

Data sheet prepared 6/19/07