

D1500.S9 / Lot No. 0610384

Beagle Dog Liver S9

Untreated, Female, Pool of 12

1.0 mL at 20 mg protein / mL

Specific content and activities^a

Content / Rate

Cytochrome P450 (nmol/mg protein)	0.130
Cytochrome b ₅ (nmol/mg protein)	0.082
7-Ethoxycoumarin <i>O</i> -dealkylation (pmol/mg protein/min) ^b	882 ± 39
Glucuronidation of 4-methylumbelliferone (nmol/mg protein/min)	151 ± 1
CDNB (nmol/mg protein/min) ^c	405 ± 8

^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: Dog; *Canis familiaris*
 Breed: Beagle
 Sex: Female
 Age: 4-9 years
 Vendor: Covance, Cumberland, VA

Dogs were laboratory animals and were housed in an AAALAC-accredited facility, which is registered as a research facility with the USDA-APHIS-AC. All animals were asymptomatic at the time of euthanasia.

Food: Nutrena (*ad libitum*)
 Water: Automatic watering system, tap water (*ad libitum*)
 Light/dark cycle: Not monitored
 Temperature: Ranges from 62°-85°F
 Humidity: Not monitored
 Cage: Indoor/outdoor run cages, plastic coated rod bottom, sanitized at least every 2 weeks



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 8/24/07

16825 West 116th St. | Lenexa KS 66219
 913.GET.P450 | fax 913.227.7100 | xenotechllc.com