

D1000.P / Lot No. 0810039

Beagle Dog Lung Microsomes
Untreated, Male, Pool of 6
0.5 mL at 10 mg protein / mL

<u>Specific content and activities</u> ^a	Content / Rate
NADPH-cytochrome <i>c</i> reductase (nmol/mg protein/min)	25.0 ± 0.3
7-Ethoxyresorufin <i>O</i> -dealkylation (pmol/mg protein/min)	9.33 ± 0.30
Glucuronidation of 4-Methylumbelliferone (nmol/mg protein/min)	< 0.1

^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.

Animal Information

Species: Dog; *Canis familiaris*
Breed: Beagle
Sex: Male
Age: > 6 months
Vendor: Covance, Cumberland, VA and Harlan

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 tissues 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 6/25/08