

## D1000.RS9 / Lot No. 0310051

Beagle Dog Renal *S9* Untreated, Male, Pool of 5 1.0 mL at 5 mg protein / mL



Specific activities <sup>a</sup>	Rate
NADPH-Cytochrome <i>c</i> Reductase (nmol/mg protein/min)	$4.65 \pm 1.05$
Lauric Acid 12-hydroxylation (pmol/mg protein/min)	$358 \pm 8$
Glucuronidation of 4-Methylumbelliferone (nmol/mg protein/min)	<1.0

<sup>&</sup>lt;sup>a</sup> Values for enzyme activities are mean + standard deviation of three or more determinations.

Subcellular fractions were prepared from whole kidney homogenate.

## **Animal Information**

Species: Dog; Canis familiaris

Breed: Beagle
Sex: Male
Age: > 6 months

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 tissue 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.

DATASHEET PREPARED 9/7/05

16825 West 116th Street, Lenexa, Kansas 66219 Phone: 913-GET-P450 www.xenotechllc.com Fax: 913-227-7100