

D1000.RS9 Lot No. 1010362

Beagle Dog Kidney S9

Untreated, Male, Pool of 5

1.0 mL at 5 mg protein / mL

Suspension medium: 50 mM Tris·HCl,
150 mM KCl,
2mM EDTA

Specific activities ^a	Rate
NADPH-cytochrome <i>c</i> reductase (nmol/mg protein/min)	11.0 ± 0.3
Lauric Acid 12-hydroxylation (pmol/mg protein/min)	549 ± 14
4-Methylumbelliferone glucuronidation (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 were determined at a single substrate concentration and are mean ± standard deviation of three or more determinations.

Animal Information

Species: Dog
Strain: Beagle
Sex: Male
Age: > 6 months
Vendor: Covance, Cumberland, VA

Animals were housed in an AAALAC-accredited facility and allowed to acclimate ≥ seven days before use.

Food: Nutrena (*ad libitum*)
Water: Automatic watering system, tap water (*ad libitum*)
Light/dark cycle: Not monitored
Temperature: Ranges from 62°-82°F
Humidity: Not monitored
Cage: Indoor/outdoor run cages, plastic coated rod bottom, sanitized at least every 2 weeks

**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 December 2010