

## D1000 Lot No. 2110225

Beagle Dog Liver Microsomes  
 Untreated, Male, Pool of 9  
 0.5 mL at 20 mg protein / mL  
 Suspension medium: 250 mM sucrose

Specific Content and Enzyme Activities		Content / Rate
Cytochrome P450 content	(nmol/mg protein)	0.587
Cytochrome b <sub>5</sub> content	(nmol/mg protein)	0.324
NADPH-cytochrome c reductase	(nmol/mg protein/min)	65.4 ± 6.6
7-Ethoxycoumarin O-dealkylation	(pmol/mg protein/min)	1990 ± 120

Values for enzyme activities were determined at a single substrate concentration and are mean ± standard deviation of three or more determinations.

To measure cytochrome P450 (CYP) activity, liver microsomes (50 µg/mL) were incubated in triplicate at 37 ± 2°C for 10 minutes in potassium phosphate buffer (50 mM, pH 7.4), containing MgCl<sub>2</sub> (3.0 mM), EDTA (1.0 mM), NADP (1.0 mM), glucose-6-phosphate (5.0 mM), glucose-6-phosphate dehydrogenase (1 Unit/mL) and 7-ethoxycoumarin (500 µM), at the final concentrations indicated. Metabolite formation was determined by validated LC-MS/MS methods with deuterated metabolites as internal standards.

### Animal Information

Species: Dog; *Canis familiaris*  
 Strain: Beagle  
 Sex: Male  
 Age: >6 months  
 Vendor: Envigo, 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. These data are not to be reproduced, published or distributed without the express written consent of XenoTech.

Datasheet prepared 12 August 2021