

## D1000 Lot No. 2310325

Beagle Dog Liver Microsomes Untreated, Male, Pool of 5 0.5 mL at 20 mg protein / mL

Suspension medium: 250 mM sucrose



Specific Content and Enzyme Activities		Content / Rate
Cytochrome P450 content Cytochrome b <sub>5</sub> content	(nmol/mg protein) (nmol/mg protein)	0.621 0.289
NADPH-cytochrome <i>c</i> reductase 7-Ethoxycoumarin <i>O</i> -dealkylation	(nmol/mg protein/min) (pmol/mg protein/min)	96.5 ± 0.6 2700 ± 230

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 ( $50 \,\mu\text{g/mL}$ ) were incubated in triplicate at  $37 \pm 2^{\circ}\text{C}$  for 10 minutes in potassium phosphate buffer ( $50 \, \text{mM}$ , pH 7.4), containing MgCl<sub>2</sub> ( $3.0 \, \text{mM}$ ), EDTA ( $1.0 \, \text{mM}$ ), NADP ( $1.0 \, \text{mM}$ ), glucose-6-phosphate ( $5.0 \, \text{mM}$ ), glucose-6-phosphate ( $5.0 \, \text{mM}$ ), glucose-6-phosphate ( $5.0 \, \text{mM}$ ), and 7-ethoxycoumarin ( $5.00 \, \mu\text{M}$ ), 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: Marshall, North Rose, NY

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 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.