

**D1073 / Lot No. 080100D**

Beagle Dog Liver Microsomes
Saline-treated
Male, Pool of 2
0.5 mL at 20 mg protein / mL

Specific content and activities	Content / Rate
Cytochrome P450 (nmol/mg protein)	0.376
Cytochrome b ₅ (nmol/mg protein)	0.235
NADPH-cytochrome <i>c</i> reductase (nmol/mg protein/min)	42.7

Background: Liver microsomes from male dogs treated with saline are intended to serve as vehicle-treated controls for studies with liver microsomes from male dogs treated with certain P450 enzyme inducers, such as phenobarbital and clofibrac acid.

Animal Information

Species:	Dog	Treatment:	Saline (0.9% NaCl irrigation solution)
Strain:	Beagle	Regimen:	5 mL/kg body weight once per day on days 1-4, liver harvested and snap-frozen on day 5
Sex:	Male		
Age:	Sexually mature		

**Store at -80°C**

CAUTION: Although strict measures are taken to ensure that livers 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.

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 expressed written consent of XENOTECH, LLC.