



D1098 / Lot No. 072800A
Beagle Dog Liver Microsomes
Corn oil-treated
Male, Pool of 2
0.5 mL at 20 mg protein / mL

Specific content and activities	Content / Rate
Cytochrome P450 (nmol/mg protein)	0.430
Cytochrome b ₅ (nmol/mg protein)	0.211
NADPH-cytochrome <i>c</i> reductase (nmol/mg protein/min)	37.8

Background: Liver microsomes from male dogs treated with corn oil are intended to serve as vehicle-treated controls for studies with liver microsomes from male dogs treated with certain P450 enzyme inducers, such as β -naphthoflavone, and rifampin.

Animal Information

Species:	Dog	Treatment:	Corn Oil
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.