

**D1095 / Lot No. 0510078**

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

Specific content and activities	Content / Rate
Cytochrome P450 (nmol/mg protein)	0.865
Cytochrome b <sub>5</sub> (nmol/mg protein)	0.265
NADPH-cytochrome <i>c</i> reductase (nmol/mg protein/min)	61.4
Testosterone 6 $\beta$ -hydroxylation (pmol/mg protein/min)	1370 <sup>a</sup>

<sup>a</sup> Fold induction: ~2.5 -fold increase over control microsomes

**Background:** Treatment of male dogs with rifampin causes a marked induction (>2-fold) of liver microsomal CYP3A levels, which is associated with an increase in testosterone 6 $\beta$ -hydroxylation. Liver microsomes from corn oil-treated dogs were used as a control.

<u>Animal Information</u>			
Species:	Dog	Treatment:	Rifampin
Strain:	Beagle	Vehicle:	Corn Oil
Sex:	Male	Regimen:	10 mg/kg body weight once per day on days 1-4, liver was harvested and snap-frozen on day 5
Age:	Sexually mature		

**Store at -80°C***For in vitro use only*

**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.

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.

DATA SHEET PREPARED 6/28/05