

R1078 Lot No. 2010127

Sprague Dawley (SD) Rat Liver Microsomes
 Phenobarbital-treated
 Male, Pool of 12
 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)	1.258
Cytochrome b ₅ content	(nmol/mg protein)	0.623
NADPH-cytochrome <i>c</i> reductase	(nmol/mg protein/min)	188 ± 2
7-Pentoxoresorufin O-dealkylation	(pmol/mg protein/min)	2120 ± 100

Background: Treatment of rats with the peroxisome proliferator, phenobarbital, causes a marked induction (>10-fold) of liver microsomal CYP2B levels, which is associated with an increase in 7-pentoxoresorufin O-dealkylation.

Animal Information

Species:	Rat	Treatment:	Phenobarbital
Strain:	* IGS Sprague Dawley	Source:	Sigma (Cat. No. P-5178)
Sex:	Male	Vehicle:	Saline, pH adjusted to 7 with NaOH
Age:	~ 8 weeks	Concentration:	16 mg/mL
Vendor:	Charles River, Raleigh, NC	Regimen:	80 mg/kg body weight once per day on days 1-4, liver microsomes prepared on day 5

*International Genetic Standard

Animals were housed in an AAALAC-accredited facility and allowed to acclimate ≥ seven days before use.

Food:	Purina 5L79 (<i>ad libitum</i>)
Water:	Automatic watering system (<i>ad libitum</i>)
Light/dark cycle:	5:00 am - 5:00 pm, light; 5:00 pm - 5:00 am, dark (12-hour light/dark)
Temperature:	70°F ± 2°F
Humidity:	30-70 %
Bedding:	Beta Chip (hardwood), NEPCO, Warrensburg, NY
Cage:	Polycarbonate Shoebox Cage, conventional cage



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 02 November 2020