

R1081.S9 Lot No. 1510257

Sprague Dawley (SD) Rat Liver S9
 β-Naphthoflavone & Phenobarbital-treated

Male, Pool of 50

1.0 mL at 20 mg protein / mL

Suspension medium: 50 mM Tris·HCl, 150 mM KCl, 2 mM EDTA

<i>Specific Content</i>		<i>Content</i>
Cytochrome P450 content	(nmol/mg protein)	0.539
Cytochrome b ₅ content	(nmol/mg protein)	0.156

Background: Dual treatment of male rats with β-naphthoflavone and the peroxisome proliferator, phenobarbital, causes a marked induction (>10-fold) of liver microsomal CYP1A and CYP2B enzymes. This dual treatment produces a similar response as Aroclor 1254, a known mixed inducer. Liver microsomes from corn oil-treated rats (Cat. No. R1098) and saline-treated rats (Cat. No. R1073) were used as controls. The results confirm the anticipated induction of both CYP1A and CYP2B activity.

Animal Information

Species:	Rat	Treatment:	β-Naphthoflavone (BNF) and Phenobarbital (PB)
Strain:	* IGS Sprague Dawley	Source:	Sigma (BNF and PB)
Sex:	Male	Vehicle:	Corn oil (BNF) and Saline (PB)
Age:	~ 8 weeks	Concentration:	BNF = 20 mg/mL, PB = 16 mg/mL
Vendor:	Charles River, Raleigh, NC	Regimen:	BNF = 100 mg/kg body weight single injection on days 1-4, livers harvested on day 5 PB = 80 mg/kg body weight single injection on days 1-4, livers harvested 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 14 December 2015