

R1083 Lot No. 1110154

Sprague Dawley (SD) Rat Liver Microsomes

β -naphthoflavone-treated, Male, Pool of 25

0.5 mL at 20 mg protein / mL

Suspension medium: 250 mM sucrose

Specific content and activities ^a

Content / Rate

Cytochrome P450	(nmol/mg protein)	1.292
Cytochrome b ₅	(nmol/mg protein)	0.678
NADPH-cytochrome <i>c</i> reductase	(nmol/mg protein/min)	148 ± 6
7-Ethoxyresorufin <i>O</i> -dealkylation	(pmol/mg protein/min)	3710 ± 300

Background: Treatment of male rats with β -naphthoflavone causes a marked induction (>10-fold) of liver microsomal CYP1A levels, which is associated with an increase in 7-ethoxyresorufin *O*-dealkylation. Liver microsomes from corn oil-treated rats (Cat. No. R1098) were used as a control. The results confirm the anticipated induction of CYP1A activity.

^a Characterization is performed when the first lot of a product from a given subcellular fraction (e.g., S9) is prepared. Subsequent lots are subject to a verification test only. Values for enzyme activities were determined at a single substrate concentration and are mean ± standard deviation of three or more determinations.

Animal Information

Species:	Rat	Treatment:	β -naphthoflavone
Strain:	* IGS, Sprague Dawley	Source:	Sigma (Cat. No. N-3633)
Sex:	Male	Vehicle:	Corn Oil
Age:	~8 weeks	Concentration:	20 mg/mL
Vendor:	Charles River, Raleigh, NC	Regimen:	100 mg/kg body weight once per day on days 1-4, livers harvested on day 5

* International Genetic Standard

Animals were housed in an AAALAC-accredited facility and allowed to acclimate \geq 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, LLC. These data are not to be reproduced, published or distributed without the express written consent of XENOTECH, LLC.

Datasheet prepared 28 September 11