

## R1078 Lot No. 1010045

Sprague Dawley (SD) Rat Liver Microsomes

Phenobarbital-treated, Male, Pool of 50

0.5 mL at 20 mg protein / mL

Suspension medium: 250 mM sucrose

### Specific content and activities <sup>a</sup>

### Content / Rate

Cytochrome P450	(nmol/mg protein)	1.85
Cytochrome b <sub>5</sub>	(nmol/mg protein)	0.629
NADPH-cytochrome <i>c</i> reductase	(nmol/mg protein/min)	202 ± 16
7-Pentoxoresorufin <i>O</i> -dealkylation (pmol/mg protein/min)		2280 ± 20

**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.

<sup>a</sup> 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:	Phenobarbital
Strain:	* IGS, Sprague Dawley	Source:	Sigma Chemical Co. (Cat. No. P-5178)
Sex:	Male	Vehicle:	Saline
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:	Harlan Teklad Rodent Chow #8604 ( <i>ad libitum</i> )
Water:	Automatic watering system ( <i>ad libitum</i> )
Light/dark cycle:	6:00 am - 6:00 pm, light; 6:00 pm - 6:00 am, dark (12-hour light/dark)
Temperature:	72°F ± 3°F
Humidity:	45-55 %
Bedding:	Cell-Sorb Plus (gypsum treated paper product), A&W Products, New Philadelphia, OH
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 16 FEBRUARY 11