

R1088 Lot No. 1110155

Sprague Dawley (SD) Rat Liver Microsomes

Isoniazid-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)	0.554
Cytochrome b ₅	(nmol/mg protein)	0.768
NADPH-cytochrome <i>c</i> reductase	(nmol/mg protein/min)	173 ± 37
4-Nitrophenol hydroxylation	(pmol/mg protein/min)	2840 ± 170

Background: Treatment of male rats with Isoniazid causes a 3- to 5-fold induction of liver microsomal CYP2E levels, which is associated with an increase in 4-nitrophenol hydroxylation, although P450 enzymes other than CYP2E1 can catalyze these same reactions.

^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:	Isoniazid
Strain:	* IGS, Sprague Dawley	Source:	Sigma (Cat. No. I-3377)
Sex:	Male	Vehicle:	Saline
Age:	~8 weeks	Concentration:	40 mg/mL
Vendor:	Charles River, Raleigh, NC	Regimen:	200 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 ≥ 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 05 October 11