



R1088 / Lot No. 0510064
 Sprague Dawley Rat Liver Microsomes
 Isoniazid-treated, Male, Pool of 25
 0.5 mL at 20 mg protein / mL

Specific content and activities
Content / Rate

Cytochrome P450 (nmol/mg protein)	1.01
Cytochrome b ₅ (nmol/mg protein)	0.791
NADPH-cytochrome <i>c</i> reductase (nmol/mg protein/min)	227 ± 10
4-nitrophenol hydroxylation (nmol/mg protein/min)	3.13 ± 0.08

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 and chlorzoxazone 6-hydroxylation, although P450 enzymes other than CYP2E1 can catalyze these same reactions.

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, liver microsomes prepared on day 5

* International Genetic Standard

Rats were laboratory animals and were housed in an AAALAC-accredited facility, which is registered as a research facility with the USDA-APHIS-AC. They were 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

For in vitro use only

CAUTION: Although strict measures are taken to ensure that livers obtained from laboratory animals do not harbor infectious diseases, we recommend that all animal products be handled as potential biohazards and universal precautions be followed.

These data were generated by and are the property of XENOTECH, LLC. These data are not to be reproduced, published or distributed without the expressed written consent of XENOTECH, LLC.

DATA SHEET PREPARED 8/24/05