



# **R1098 / Lot No. 0510066**

Sprague Dawley Rat Liver Microsomes  
Corn Oil-treated, Male, Pool of 25  
0.5 mL at 20 mg protein / mL

Specific content and activities	Content / Rate
Cytochrome P450 (nmol/mg protein)	0.802
Cytochrome b <sub>5</sub> (nmol/mg protein)	0.498
NADPH-cytochrome <i>c</i> reductase (nmol/mg protein/min)	184 ± 15

**Background:** Liver microsomes from male rats treated with corn oil are intended to serve as vehicle-treated controls for studies with liver microsomes from male rats treated with certain P450 enzyme inducers, such as  $\beta$ -naphthoflavone, dexamethasone, 3-methylcholanthrene, Aroclor 1254, pregnenolone-16 $\alpha$ -carbonitrile and perfluorodecanoic acid.

Note: XenoTech also offers liver microsomes from untreated rats (R1000) and rats treated with saline (R1073).

## **Animal Information**

Species:	Rat	Treatment:	Corn Oil
Strain:	Sprague Dawley	Source:	Sigma (Cat. No. C-8267)
Sex:	Male	Volume injected:	5 mL/kg body weight
Age:	~8 weeks	Regimen:	Once per day on days 1-4, liver microsomes prepared on day 5
Vendor:	Charles River, Raleigh, NC		

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  $\geq$  seven days before use.

Food:	Purina Rodent Laboratory Chow #5001 ( <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 $\pm$ 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 7/14/05