

**P2573 Lot No. PR10021**

Cynomolgus Monkey Liver Microsomes

Saline-treated, Female, Pool of 2

0.5 mL at 10 mg protein / mL

Suspension medium: 250 mM sucrose

Specific content and activities <sup>a</sup>		Content / Rate
Cytochrome P450	(nmol/mg protein)	0.847
Cytochrome b <sub>5</sub>	(nmol/mg protein)	0.259
NADPH-cytochrome <i>c</i> reductase	(nmol/mg protein/min)	101 ± 13

**Background:** Liver microsomes from female monkeys treated with saline are intended to serve as vehicle-treated controls for studies with liver microsomes from female monkeys treated with certain P450 enzyme inducers, such as phenobarbital, rifampin and omeprazole.

**Animal Information**

Species:	Monkey	Treatment:	Saline (0.9% NaCl irrigation solution)
Strain:	Cynomolgus	Supplier:	Baxter (Cat. No. 2F7124)
Sex:	Female	Regimen:	Treatment once per day on days 1-4,
Age:	Sexually mature		tissue harvested on day 5.
Vendor:	WIL Research		

Monkeys were housed in an AAALAC-accredited facility

Imported animals were quarantined for one month prior to shipment into the United States to reduce the risk of importing Ebola virus-infected monkeys. All animals were under veterinary care and were asymptomatic at the time of euthanasia. All of the monkeys tested negative for Simian Retrovirus. None of the animals examined tested positive for any other infectious agents.

**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 10 August 2010