



EAGLE
BIOSCIENCES

Monoclonal Anti-Rat Cytochrome P450 (2B1)

Product Number: CYT00-A100

Aliquot: 100 μ L

Storage: - 20°C

PRODUCT SPECIFICATION SHEET

Type:	Monoclonal Mouse IgG
Specificity:	Identification of 2B1 by immunoassay or Western blots.
Applications:	Immunoassay or Western blots
Contents:	100 μ L per vial. Contains sufficient purified monoclonal mouse IgG for at least 10 Western blots.
Working Dilution:	A dilution of 1/1000 for Western blots will yield a strong band with 10 μ g of rat microsomes using Alkaline Phosphatase.
Storage:	Stable for several weeks at 4°, several months at -20°, and indefinitely at -80°. Do not freeze and thaw repeatedly. Refreezing the product may result in diminishment or loss of activity.

Product Developed and Manufactured in the USA

EAGLE BIOSCIENCES, INC.
20A NW Blvd., Suite 112
Nashua, NH 03063

Phone: 617-419-2019
Fax: 617-419-1110
Email: info@EagleBio.com
www.EagleBio.com