



**EAGLE**  
BIOSCIENCES

## Monoclonal Anti-Rat Cytochrome P450 (2B1, 2B2)

Product Number: CYT00-A1000

Aliquot: 1 mL

Storage: -20°C

### PRODUCT SPECIFICATION SHEET

**Immunoglobulin**

**Type:** [Monoclonal Mouse IgG.](#)

**Specificity:** Identification of 2B1 and 2B2 by immunoassay or Western blots.

**Applications:** Immunoassay or Western blots.

**Contents:** 1mL per vial. Contains sufficient purified monoclonal mouse IgG for at least 10 Western blots.

**Working Dilution:** A dilution of 1/100 for Western blots will yield a strong band with 10 µg of Phenobarbital induced rat microsomes using Alkaline Phosphatase.

**Storage:** Stable for several weeks at 4°C, several months at -20°C, and indefinitely at -80°C. 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](mailto:info@EagleBio.com)  
[www.EagleBio.com](http://www.EagleBio.com)