



**EAGLE**  
BIOSCIENCES

## Anti-Rabbit Cytochrome P450 1A1/1A2

Product Number: CPA00-A1000

Aliquot: 1 mL

Storage: -20°C

### PRODUCT SPECIFICATION SHEET

#### Immunoglobulin

**Type:**

Polyclonal goat IgG.

**Immunogen:**

Purified rabbit CYP450 1A1/1A2.

**Specificity:**

Reacts with both forms in liver microsomes prepared from rabbits or rats that were induced to express multiple forms of P450's.

**Applications:**

Immunoassay, Immunoprecipitation or Western Blots.

**Contents:**

1 ml of goat antiserum with 0.05% sodium azide.

**Working Dilution:**

A dilution of 1/100 for Western blots will yield a positive result with 0.2 µg of recombinant human cytochrome P450 1A1 and/ or 1A2 using horseradish conjugated rabbit anti-goat antibodies according to the manufacturer's instructions.

**Storage:**

Stable at 4°C for 6 to 8 weeks, several months at -20°C, and indefinitely at -80°C. Upon thawing it is best to aliquot for individual usage. Repeated freezing and thawing may result in diminishment or loss of activity.

**References:**

1. Guengerich, F.P. *et al.* (1982) *Biochemistry* **21**: 6019
2. Bellward, G.D., *et al.* (1988) *Mol. Pharmacol.* **33**: 140

*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)