



EAGLE
BIOSCIENCES

Anti 2B4 Cytochrome P450

Product Number: ACY00-A1000

Aliquot: 1 mL

Storage: - 80°C

PRODUCT SPECIFICATION SHEET

Immunoglobulin Type:	Polyclonal goat IgG.
Specificity:	Reacts as a single band with liver microsomes prepared from rabbits that were induced to express multiple forms of P450's.
Applications:	Immunoassay or Western blots.
Contents:	1 mL of goat antiserum containing 0.05% sodium azide.
Working Dilution:	A dilution of 1/100 for Western blots will yield a positive result with 0.2 µg of purified rabbit cytochrome P450 2B4 using horseradish peroxidase conjugated rabbit anti-goat antibodies according to the manufacturer's instructions. Diluted serum can be used a minimum of three times.
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:	<ol style="list-style-type: none">1. Guengerich, F.P., <i>et al.</i> (1982) <i>Biochemistry</i> 21: 60192. Kim, S.G., <i>et al.</i> (1988) <i>J. Pharmacol. Exp. Ther.</i> 246: 11753. Bellward, G.D., <i>et al.</i> (1988) <i>Mol. Pharmacol.</i> 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
www.EagleBio.com