



EAGLE
BIOSCIENCES

Anti-Protein Kinase C Eta (η)

Product Number: KCH00-A100

Aliquot: 100 μ L

Storage: -20°C

PRODUCT SPECIFICATION SHEET

Immunoglobulin

Type: Polyclonal Rabbit IgG

Specificity: Prepared against synthetic peptide corresponding to a distinct C terminal region, from rat protein kinase C η and based on known homologies is expected to react with the PKC ortholog in most mammalian species.

Applications: Intended for use in Western blots, ELISA, Immunocytochemistry and Immunohistochemical techniques for fresh or frozen sections, paraffin embedded or fixed tissues have not been tested.

dilutions: For Immunohistochemistry, immunocytochemistry and Western blotting using alkaline phosphatase or horseradish peroxidase, a dilution of 1:250 is recommended. With a more sensitive detection system, such as enhanced chemiluminescence or ELISA, the antibody may be diluted further.

Contents per vial: 100 μ L of rabbit antiserum containing 0.05% sodium azide.

Storage: Stable for several months at 4°C, up to a year at -20°C, and indefinitely at -80° C. Avoid repeated freeze/thaw cycles.

References:

1. Wetsel, W. C., et al., *J. Cell Biol.* **117**: 21-133 (1992)
2. *JBC*, **268** : 20110-20115

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