



**EAGLE**  
BIOSCIENCES

## Anti-Protein Kinase C Delta ( $\delta$ )

Product Number: KCD00-A100

Aliquot: 100  $\mu$ L

Storage: -20°C

### PRODUCT SPECIFICATION SHEET

#### Immunoglobulin

**Type:** Polyclonal Rabbit IgG

**Specificity:** Prepared against synthetic peptide corresponding to a distinct C terminal region (SFVNPKEYEQFLE residues 726-737) from rat protein kinase C $\delta$  and based on known homologies is expected to react with 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.

**Recommended 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  $\mu$ 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](mailto:info@EagleBio.com)  
[www.EagleBio.com](http://www.EagleBio.com)