



EAGLE
BIOSCIENCES

Anti-Protein Kinase C δ

Catalog Number: PKC00-A1000

Aliquot: 1000 μ L

Storage: -70°C

PRODUCT SPECIFICATION SHEET

DESCRIPTION:	This polyclonal antibody is produced in rabbit and is provided as antiserum specific for Protein Kinase C δ (PKC δ).
MOLECULAR WEIGHT:	77.5 kDa Calculated
CONCENTRATION:	0.68 mg/mL
ISOTYPE:	NA
STORAGE BUFFER:	Phosphate Buffered Saline; 0.05% Sodium Azide
STORAGE:	-70°C. AVOID REPEATED FREEZE THAW CYCLES. Has demonstrated stability at 4 °C for greater than 1 month.
IMMUNOGEN:	KLH conjugated synthetic peptide corresponding to amino acids 658- 676 of PKC δ . Sequence: ⁶⁵⁸ S-A-F-A-G-F-S-F-V-N-P-K-F-E-L-L-E-D ⁶⁷⁶
PURIFICATION:	NA
APPLICATIONS:	Western Blotting: 1:100 Dilution
SPECIES SPECIFICITY:	Human, mouse and rat among other mammalian species.
REFERENCE:	JBC (1993) 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