



EAGLE
BIOSCIENCES

Anti-Protein Kinase C Alpha α

Catalog Number: PKA00-A1000

Aliquot: 1000 μ L

Storage: -70°C

PRODUCT SPECIFICATION SHEET

DESCRIPTION:	This polyclonal antibody is produced in rabbit and is provided as an IgG presentation for Protein Kinase C- α (PKC α).
MOLECULAR WEIGHT:	76.8 kDa Calculated
CONCENTRATION:	1 mg/mL
ISOTYPE:	IgG
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 662- 672 of PKC α . Sequence: ⁶⁶³ Q-F-V-H-P-I-L-Q-S-S-V ⁶⁷²
PURIFICATION:	Ammonium Sulfate Fractionation
APPLICATIONS:	Western Blotting: 1-5 μ g/mL
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