
Pka-C1 Polyclonal Antibody

(Catalog # A64270)

Background

cAMP-dependent protein kinase complex, cell body, cytoplasm, dendrite, plasma membrane, cAMP-dependent protein kinase activity, protein serine/threonine kinase activity, anesthesia-resistant memory, anterior/posterior pattern specification, imaginal disc, behavioral response to ethanol

Description

Pka-C1 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Liquid. 0.03% Proclin 300. 50% Glycerol, 0.01M PBS, PH 7.4.

Specificity

Drosophila melanogaster

Isotype

IgG

Uniprot ID

P12370

Purification

Protein G purified

Immunogen

Recombinant Drosophila melanogaster cAMP-dependent protein kinase catalytic subunit 1 protein (2-353AA)

Storage

Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

cAMP-dependent protein kinase catalytic subunit, PKA C, Pka-C1, CdkA, DC0, CG4379

Application

ELISA; Recommended dilution: ELISA 1:500-1:2000
