
PRKRA Polyclonal Antibody

(Catalog #A72895)

Background

This gene encodes a protein kinase activated by double-stranded RNA which mediates the effects of interferon in response to viral infection. Mutations in this gene have been associated with dystonia. Alternative splicing results in multiple transcript variants.

Description

PRKRA Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Buffer: PBS with 0.02% sodium azide, 50% glycerol, pH7.3.

Specificity

Mouse

Isotype

IgG

Uniprot ID

O75569

Purification

Affinity Purification

Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 1-313 of human PRKRA (NP_003681.1).

Storage

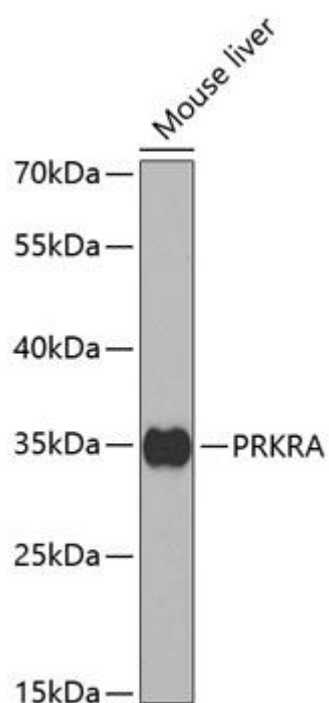
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

Alternative Names

PRKRA; DYT16; HSD14; PACT; RAX; protein activator of interferon induced protein kinase EIF2AK2

Application

WB, IHC, ELISA; Recommended dilution: WB: 1:500-1:2000, IHC: 1:50-1:200, ELISA - recommended starting concentration is 1 µg/mL. Please optimize the concentration based on your specific assay requirements.



Western blot analysis of extracts of mouse liver, using
PRKRA Polyclonal Antibody at 1:1000 dilution.
Secondary antibody: HRP Goat Anti-Rabbit IgG (H+L) at
1:10000 dilution.
Lysates/proteins: 25ug per lane.
Blocking buffer: 3% nonfat dry milk in TBST.
Exposure time: 90s.