

GRK4 Polyclonal Antibody

(Catalog # A70500)

Background

This gene encodes a member of the guanine nucleotide-binding protein (G protein)-coupled receptor kinase subfamily of the Ser/Thr protein kinase family. The protein phosphorylates the activated forms of G protein-coupled receptors thus initiating its deactivation. This gene has been linked to both genetic and acquired hypertension. Several transcript variants encoding different isoforms have been found for this gene.

Description

GRK4 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human, Mouse, Rat

Isotype

IgG

Uniprot ID

P32298

Purification

Affinity Purification

Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 359-578 of human GRK4 (NP_892027.2).

Storage

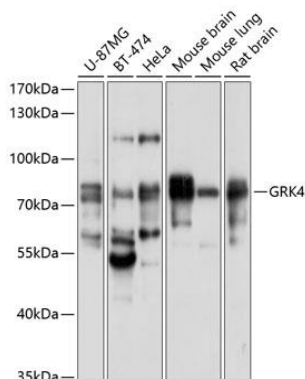
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

Alternative Names

GRK4; GPRK2L; GPRK4; GRK4a; IT11; G protein-coupled receptor kinase 4

Application

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



Western blot analysis of extracts of various cell lines, using GRK4 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: 1s.