

CHCHD3 Polyclonal Antibody

(Catalog # A73490)

Background

The protein encoded by this gene is an inner mitochondrial membrane scaffold protein. Absence of the encoded protein affects the structural integrity of mitochondrial cristae and leads to reductions in ATP production, cell growth, and oxygen consumption. This protein is part of the mitochondrial contact site and cristae organizing system (MICOS). Several transcript variants encoding different isoforms have been found for this gene.

Description

CHCHD3 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

Q9NX63

Purification

Affinity Purified

Immunogen

Recombinant protein of human CHCHD3.

Storage

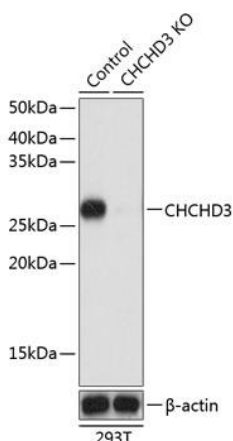
Shipped at 4°C. Store at -20°C. Avoid freeze / thaw cycles.

Alternative Names

CHCHD3; MINOS3; Mic19; PPP1R22

Application

WB, IHC; Recommended dilution: WB 1:500 - 1:2000, IHC 1:50 - 1:100



Western blot analysis of extracts from normal (control) and CHCHD3 knockout (KO) 293T cells, using CHCHD3 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: 30s.