

LAMTOR1 Polyclonal Antibody

(Catalog # A73512)

Description

LAMTOR1 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

Q6IAA8

Purification

Affinity Purified

Immunogen

Recombinant protein of human LAMTOR1.

Storage

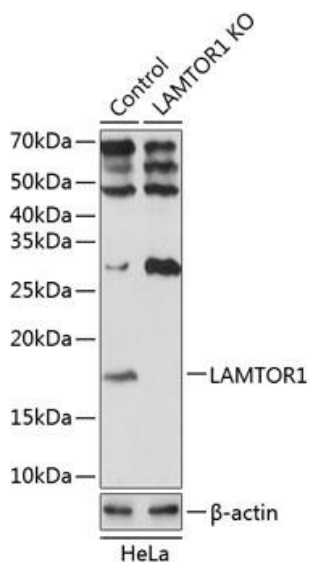
Shipped at 4°C. Store at -20°C. Avoid freeze / thaw cycles.

Alternative Names

LAMTOR1; C11orf59; PDRO; Ragulator1; p18; p27RF-Rho

Application

WB, IHC, IF; Recommended dilution: WB 1:500 - 1:1000, IHC 1:50 - 1:100, IF 1:50 - 1:100



Western blot analysis of extracts from normal (control) and LAMTOR1 knockout (KO) HeLa cells, using LAMTOR1 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.