

CLUH Polyclonal Antibody

(Catalog # A73508)

Description

CLUH Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

O75153

Purification

Affinity Purified

Immunogen

Recombinant protein of human CLUH.

Storage

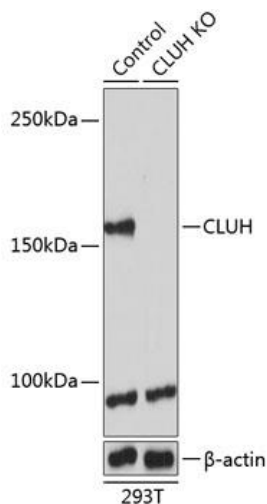
Shipped at 4°C. Store at -20°C. Avoid freeze / thaw cycles.

Alternative Names

CLUH; CLU1; KIAA0664

Application

WB, IHC, IF; Recommended dilution: WB 1:1000 - 1:4000, IHC 1:50 - 1:200, IF 1:50 - 1:200



Western blot analysis of extracts from normal (control) and CLUH knockout (KO) 293T cells, using CLUH 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.