

LYAR Polyclonal Antibody

(Catalog # A73448)

Description

LYAR 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

Q9NX58

Purification

Affinity Purified

Immunogen

Recombinant protein of human LYAR.

Storage

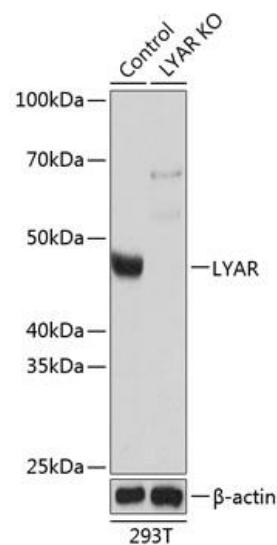
Shipped at 4°C. Store at -20°C. Avoid freeze / thaw cycles.

Alternative Names

LYAR; ZC2HC2; ZLYAR

Application

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



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