

LYAR Polyclonal Antibody

(Catalog #A72842)

Description

LYAR Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Buffer: PBS with 0.01% thimerosal, 50% glycerol, pH7.3.

Specificity

Human, Mouse, Rat

Isotype

IgG

Uniprot ID

Q9NX58

Purification

Affinity Purification

Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 10-160 of human LYAR (NP_060286.1).

Storage

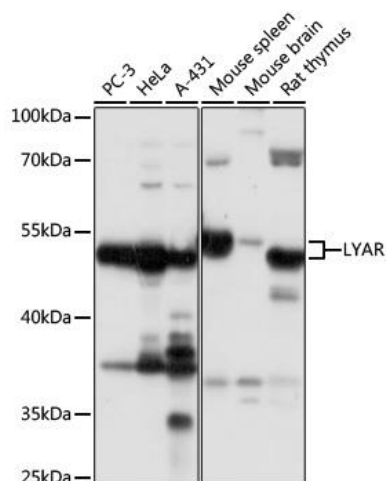
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

Alternative Names

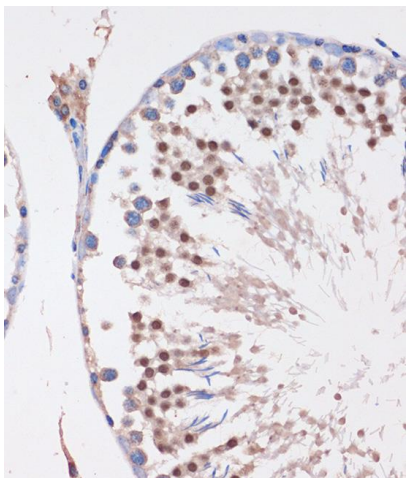
ZC2HC2, ZLYAR

Application

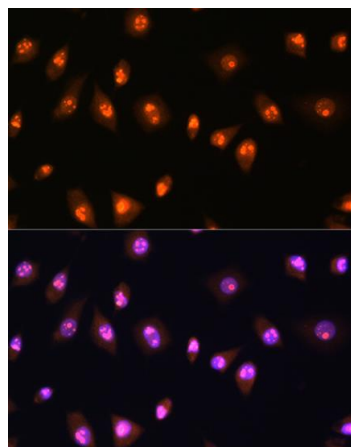
ELISA, WB, IHC, IF; Recommended dilution: WB 1:500-1:2000, IHC 1:50-1:200, IF 1:50-1:200



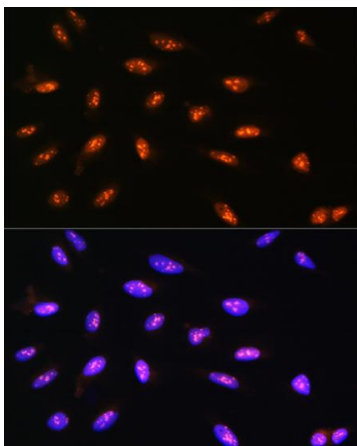
Western blot analysis of various lysates using LYAR Polyclonal Antibody at 1:1000 dilution.
Secondary antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution.
Lysates/proteins: 25µg per lane.
Blocking buffer: 3% nonfat dry milk in TBST.
Exposure time: 5s.



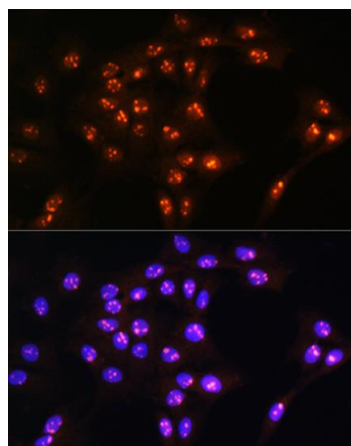
Immunohistochemistry analysis of paraffin-embedded Rat testis using LYAR Polyclonal Antibody at dilution of 1:100 (40x lens). Perform microwave antigen retrieval with 10 mM PBS buffer pH 7.2 before commencing with IHC staining protocol.



Immunofluorescence analysis of L929 cells using LYAR Polyclonal Antibody at dilution of 1:100. Secondary antibody: Cy3 Goat Anti-Rabbit IgG (H+L) at 1:500 dilution. Blue: DAPI for nuclear staining.



Immunofluorescence analysis of U2OS cells using LYAR Polyclonal Antibody at dilution of 1:100. Secondary antibody: Cy3 Goat Anti-Rabbit IgG (H+L) at 1:500 dilution. Blue: DAPI for nuclear staining.



Immunofluorescence analysis of C6 cells using LYAR Polyclonal Antibody at dilution of 1:100. Secondary antibody: Cy3 Goat Anti-Rabbit IgG (H+L) at 1:500 dilution. Blue: DAPI for nuclear staining.