
RNASE2 Polyclonal Antibody

(Catalog # A71043)

Background

The protein encoded by this gene is a non-secretory ribonuclease that belongs to the pancreatic ribonuclease family, a subset of the ribonuclease A superfamily. The protein has antimicrobial activity against viruses.

Description

RNASE2 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Mouse

Isotype

IgG

Uniprot ID

P10153

Purification

Affinity Purification

Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 28-161 of human RNASE2 (NP_002925.1).

Storage

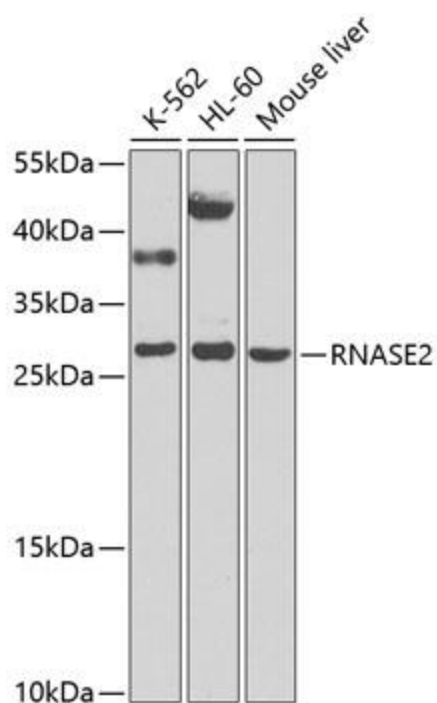
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

Alternative Names

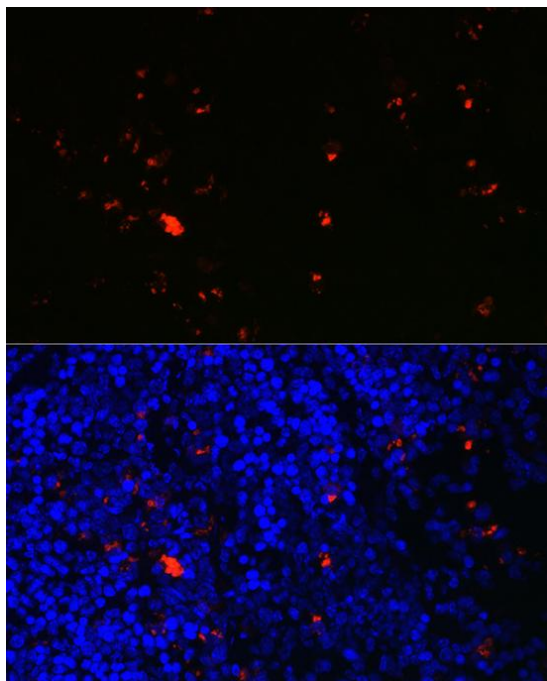
RNASE2; EDN; RAF3; RNS2; non-secretory ribonuclease

Application

WB, IF, ELISA; Recommended dilution: WB: 1:500 - 1:2000, IF: 1:50-1:200, ELISA - recommended starting concentration is 1 µg/mL. Please optimize the concentration based on your specific assay requirements.



Western blot analysis of extracts of various cell lines, using RNASE2 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: 90s.



Immunofluorescence analysis of mouse spleen using RNASE2 Polyclonal Antibody dilution of 1:100 (40x lens). Blue: DAPI for nuclear staining.