
RNASE3 Polyclonal Antibody

(Catalog #A72637)

Background

The protein encoded by this gene belongs to the pancreatic ribonuclease family, a subset of the ribonuclease A superfamily. The protein exhibits antimicrobial activity against pathogenic bacteria

Description

RNASE3 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human, Rat

Isotype

IgG

Uniprot ID

P12724

Purification

Affinity Purified

Immunogen

A synthetic peptide corresponding to a sequence within amino acids 70-150 of human RNASE3 (NP_002926.2)

Storage

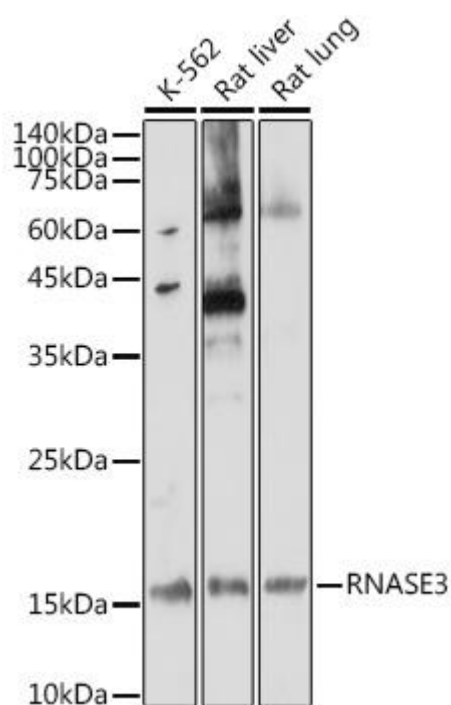
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

Alternative Names

RNASE3; ECP; RAF1; RNS3; eosinophil cationic protein

Application

WB, ELISA; Recommended dilution: WB 1:500 - 1:1000, ELISA - recommended starting concentration is 1 µg/mL. Please optimize the concentration based on your specific assay requirements.



Western blot analysis of various lysates using RNASE3 Polyclonal Antibody at 1:1000 dilution.
Secondary antibody: HRP-conjugated 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: 180s.