

RNASE2 Polyclonal Antibody

(Catalog #A52888)

Background

This is a non-secretory ribonuclease. It is a pyrimidine specific nuclease with a slight preference for U. Cytotoxin and helminthotoxin. Selectively chemotactic for dendritic cells. Possesses a wide variety of biological activities.

Description

RNASE2 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Liquid. 0.03% Proclin 300, 50% Glycerol, 0.01M PBS, PH 7.4.

Specificity

Human

Isotype

IgG

Uniprot ID

P10153

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Non-secretory ribonuclease protein (28-161AA)

Storage

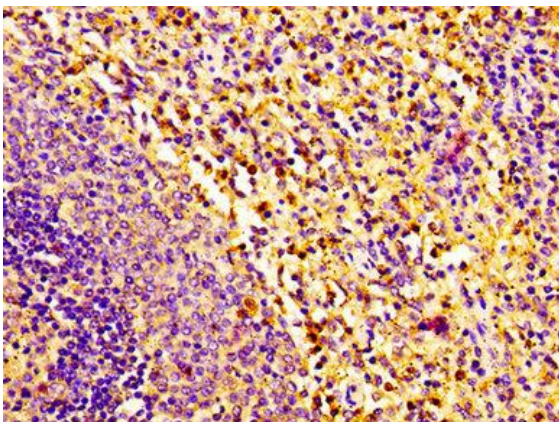
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Eosinophil-derived neurotoxin, RNase Upl-2, Ribonuclease 2, Ribonuclease US, RNASE2, EDN, RNS2

Application

ELISA, IHC; Recommended dilution: IHC:1:20-1:200



IHC image of RNASE2 Antibody diluted at 1:100 and staining in paraffin-embedded human spleen tissue performed on a Leica Bond™ system. After dewaxing and hydration, antigen retrieval was mediated by high pressure in a citrate buffer (pH 6.0). Section was blocked with 10% normal goat serum 30min at RT. Then primary antibody (1% BSA) was incubated at 4°C overnight. The primary is detected by a biotinylated secondary antibody and visualized using an HRP conjugated SP system.