
COLEC10 Polyclonal Antibody

(Catalog # A56980)

Background

Lectin that binds to various sugars: galactose > mannose = fucose > N-acetylglucosamine > N-acetylgalactosamine.

Description

COLEC10 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human

Isotype

IgG

Uniprot ID

Q9Y6Z7

Purification

Protein G purified

Immunogen

Recombinant Human Collectin-10 protein (28-277AA)

Storage

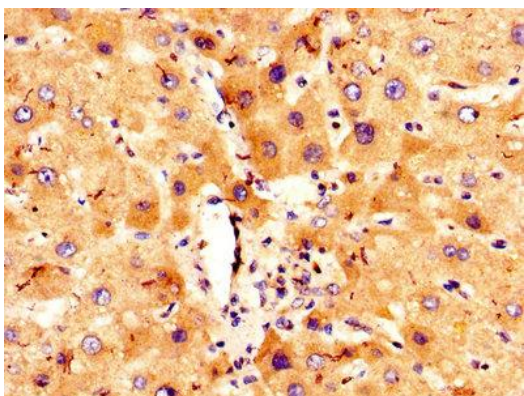
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

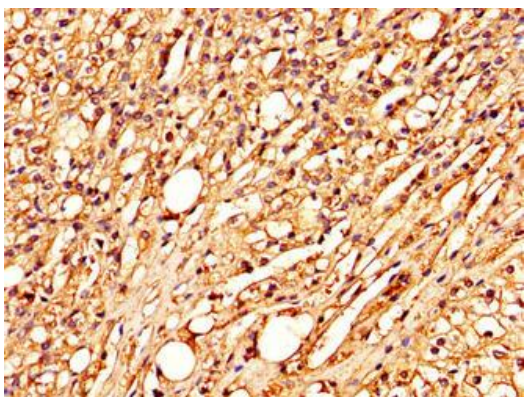
Collectin liver protein 1, CL-L1, Collectin-34, CL-34, UNQ366/PRO702, CLL1, COLEC10

Application

ELISA, IHC; Recommended dilution: IHC 1:200 - 1:500



IHC image of COLEC10 Antibody diluted at 1:300 and staining in paraffin-embedded human liver 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.



IHC image of COLEC10 Antibody diluted at 1:300 and staining in paraffin-embedded human adrenal gland 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.