
RNF138 Polyclonal Antibody

(Catalog #A60626)

Background

E3 ubiquitin-protein ligase. Together with NLK, involved in the ubiquitination and degradation of TCF/LEF. Also exhibits auto-ubiquitination activity in combination with UBE2K. May act as a negative regulator in the Wnt/beta-catenin-mediated signaling pathway.

Description

RNF138 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

Q8WVD3

Purification

>95%, Protein G purified

Immunogen

Recombinant Human E3 ubiquitin-protein ligase RNF138 protein (2-245AA)

Storage

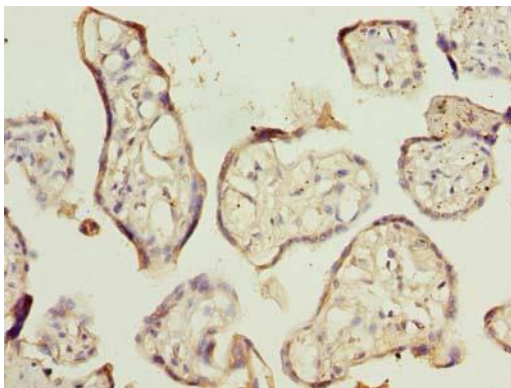
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

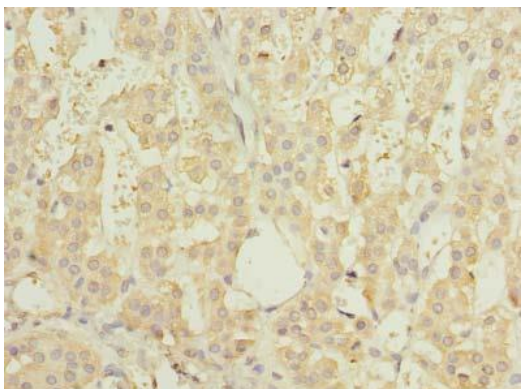
E3 ubiquitin-protein ligase RNF138 (EC:6.3.2.-), RNF138, NARF, HSD-4, HSD4, Nemo-like kinase-associated RING finger protein, NLK-associated RING finger protein, hNARF, RING finger protein 138

Application

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



Immunohistochemistry of paraffin-embedded human placenta tissue using RNF138 Polyclonal Antibody at dilution of 1:100



Immunohistochemistry of paraffin-embedded human adrenal gland tissue using RNF138 Polyclonal Antibody at dilution of 1:100