

COMMD2 Polyclonal Antibody

(Catalog # A65926)

Background

May modulate the activity of cullin-RING E3 ubiquitin ligase (CRL) complexes (PubMed:21778237). May down-regulate activation of NF-kappa-B (PubMed:15799966).

Description

COMMD2 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

Q86X83

Purification

>95%, Protein G purified

Immunogen

Recombinant Human COMM domain-containing protein 2 protein (5-199AA)

Storage

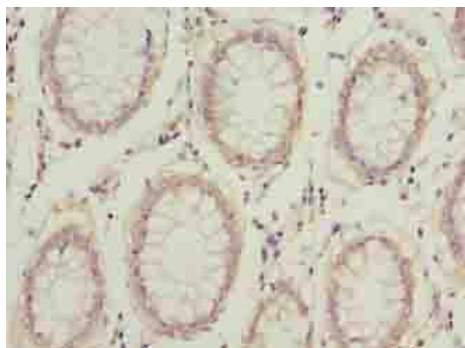
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

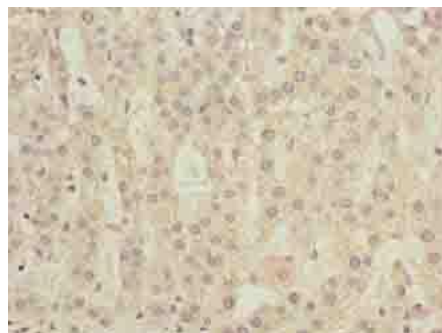
COMM domain-containing protein 2, COMMD2, HSPC042, My004

Application

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



Immunohistochemistry of paraffin-embedded human colon cancer using COMMD2 Polyclonal Antibody at dilution of 1:100



Immunohistochemistry of paraffin-embedded human adrenal gland tissue using COMMD2 Polyclonal Antibody at dilution of 1:100