

COMMD6 Polyclonal Antibody

(Catalog # A65854)

Background

May modulate activity of cullin-RING E3 ubiquitin ligase (CRL) complexes (PubMed:21778237). Down-regulates activation of NF-kappa-B. Inhibits TNF-induced NFKB1 activation.

Description

COMMD6 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

Q7Z4G1

Purification

>95%, Protein G purified

Immunogen

Recombinant Human COMM domain-containing protein 6 protein (1-85AA)

Storage

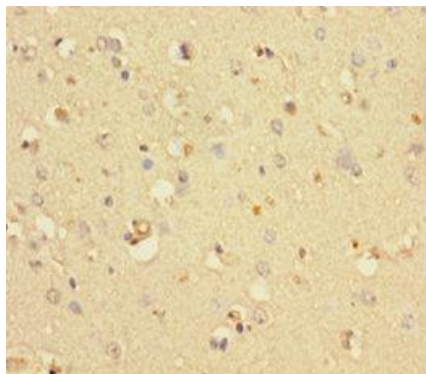
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

COMM domain-containing protein 6, COMMD6, MSTP076

Application

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



Immunohistochemistry of paraffin-embedded human brain tissue using COMMD6 Polyclonal Antibody at dilution 1:100