
CCDC79 Polyclonal Antibody

(Catalog # A66382)

Background

Meiosis-specific telomere-associated protein involved in meiotic telomere clustering, also named bouquet formation, the telomere-mediated chromosomal movement along the nuclear envelope which is crucial for homologous pairing and synapsis. Probably acts as a hub for the assembly of a conserved meiotic telomere complex required for chromosome movements via interaction with TERF1. Promotes telomere association with the nuclear envelope and deposition of the SUN-KASH/LINC complex. Also recruits cohesin to telomeres to develop structural rigidity.

Description

CCDC79 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

Q8NA31

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Telomere repeats-binding bouquet formation protein 1 protein (301-550AA)

Storage

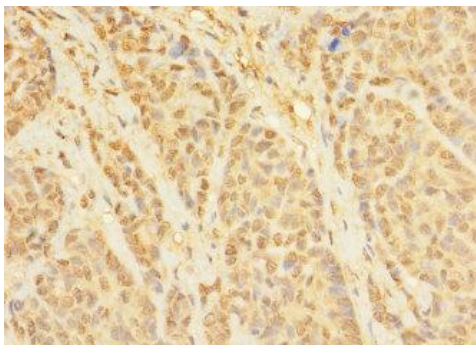
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

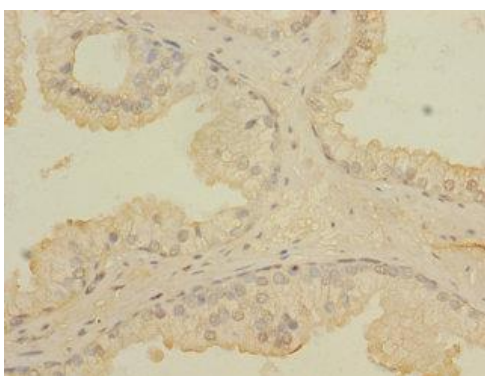
Telomere repeats-binding bouquet formation protein 1, Coiled-coil domain-containing protein 79, CCDC79, TERB1

Application

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



Immunohistochemistry of paraffin-embedded human ovarian cancer CCDC79 Polyclonal Antibody at dilution 1:100



Immunohistochemistry of paraffin-embedded human prostate cancer CCDC79 Polyclonal Antibody at dilution 1:100