

## CCDC6 Polyclonal Antibody

(Catalog # A73449)

### Background

This gene encodes a coiled-coil domain-containing protein. The encoded protein is ubiquitously expressed and may function as a tumor suppressor. A chromosomal rearrangement resulting in the expression of a fusion gene containing a portion of this gene and the intracellular kinase-encoding domain of the ret proto-oncogene is the cause of thyroid papillary carcinoma.

### Description

CCDC6 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

### Formulation

Buffer: PBS with 0.02% sodium azide, 50% glycerol, pH7.3.

### Specificity

Human, Mouse, Rat

### Isotype

IgG

### Uniprot ID

Q16204

### Purification

Affinity Purified

### Immunogen

Recombinant protein of human CCDC6.

### Storage

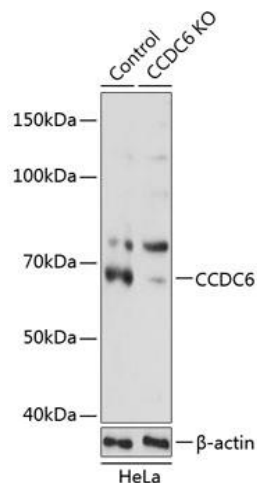
Shipped at 4°C. Store at -20°C. Avoid freeze / thaw cycles.

### Alternative Names

CCDC6; D10S170; H4; PTC; TPC; TST1

### Application

WB, IF; Recommended dilution: WB 1:500 - 1:2000, IF 1:50 - 1:200



Western blot analysis of extracts from normal (control) and CCDC6 knockout (KO) HeLa cells, using CCDC6 Polyclonal Antibody at 1:1000 dilution.  
Secondary Antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution.  
Lysates/proteins: 25ug per lane.  
Blocking buffer: 3% nonfat dry milk in TBST. Exposure time: 1s.