

## NFKBID Polyclonal Antibody

(Catalog #A66786)

### Background

May regulate the expression of IL-2, IL-6, and other cytokines through regulation on NF-kappa-B activity. Functions in the regulation of inflammatory responses. May also regulate TCR-induced negative selection of thymocytes.

### Description

NFKBID Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

### Formulation

0.03% Proclin 300, 50% Glycerol, 0.01M PBS, PH 7.4.

### Specificity

Human

### Isotype

IgG

### Uniprot ID

Q8NI38

### Purification

Protein G purified

### Immunogen

Recombinant Human NF-kappa-B inhibitor delta protein (1-150AA)

### Storage

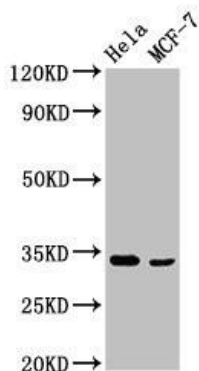
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

### Alternative Names

NF-kappa-B inhibitor delta, I-kappa-B-delta, Ikb-delta, IkappaBdelta, IkappaBNS, T-cell activation NFKB-like protein, TA-NFKBH, NFKBID, IKBNS

### Application

ELISA, WB; Recommended dilution: WB:1:500-1:2000



#### Western Blot

Positive WB detected in: HeLa whole cell lysate, MCF-7 whole cell lysate

All lanes: NFKBID Polyclonal Antibody at 3.4ug/ml

Secondary

Goat polyclonal to rabbit IgG at 1/50000 dilution

Predicted band size: 34, 51, 16 KDa

Observed band size: 34 KDa