

---

## STEAP1 Polyclonal Antibody

(Catalog # A54181)

---

### Background

E3 ubiquitin-protein ligase that mediates ubiquitination of Delta receptors, which act as ligands of Notch proteins. Positively regulates the Delta-mediated Notch signaling by ubiquitinating the intracellular domain of Delta, leading to endocytosis of Delta receptors. Probably mediates ubiquitination and subsequent proteasomal degradation of DAPK1, thereby antagonizing anti-apoptotic effects of DAPK1 to promote TNF-induced apoptosis. By similarity. Mediates 'Lys-63'-linked polyubiquitination of TBK1, which probably participates in kinase activation.

### Description

STEAP1 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

Q9UHE8

### Purification

>95%, Protein G purified

### Immunogen

Recombinant Human Metalloreductase STEAP1 protein (3-69AA)

### Storage

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

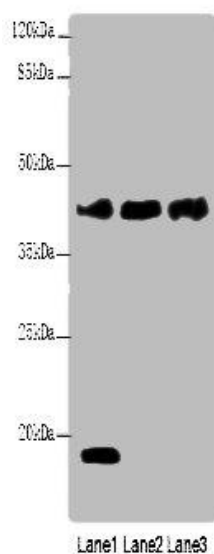
### Alternative Names

DAPK-interacting protein 1, DIP-1, Mind bomb homolog 1, Zinc finger ZZ type with ankyrin repeat domain protein 2, DIP1, KIAA1323, ZZANK2, MIB1

### Application

ELISA, WB, IHC, IF; Recommended dilution: WB:1:500-1:2000, IHC:1:20-1:200, IF:1:50-1:200

---



#### Western blot

All lanes: STEAP1 Polyclonal Antibody at 1ug/ml

Lane 1: U87 whole cell lysate

Lane 2: A431 whole cell lysate

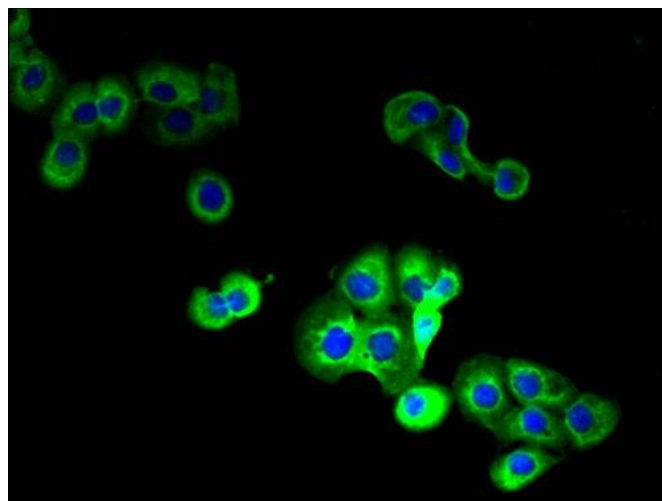
Lane 3: A549 whole cell lysate

Secondary

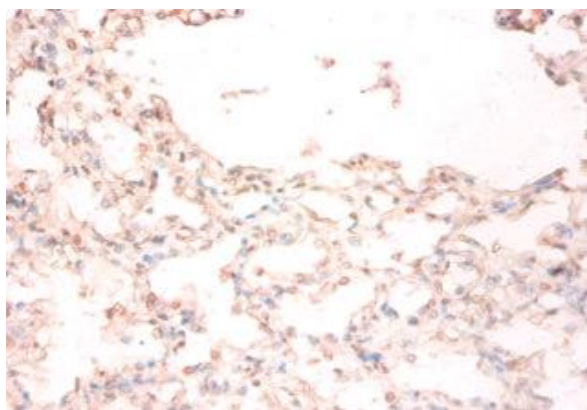
Goat polyclonal to rabbit IgG at 1/10000 dilution

Predicted band size: 40 kDa

Observed band size: 40, 18 kDa



Immunofluorescent analysis of A431 cells at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)



Immunohistochemistry of paraffin-embedded human lung at dilution of 1:50