

## BMPR1A Polyclonal Antibody

(Catalog # A50089)

### Background

On ligand binding, forms a receptor complex consisting of two type II and two type I transmembrane serine/threonine kinases. Type II receptors phosphorylate and activate type I receptors which autophosphorylate, then bind and activate SMAD transcriptional regulators. Receptor for BMP-2 and BMP-4. Positively regulates chondrocyte differentiation through GDF5 interaction.

### Description

BMPR1A 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

P36894

### Purification

Protein G purified

### Immunogen

Recombinant Human Bone morphogenetic protein receptor type-1A protein (177-532AA)

### Storage

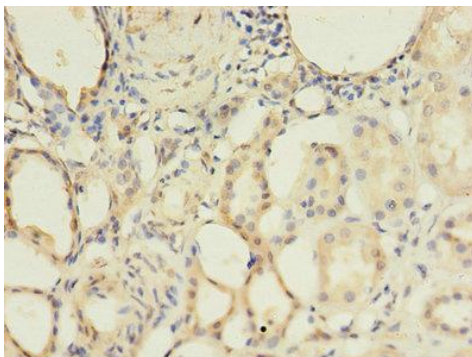
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

### Alternative Names

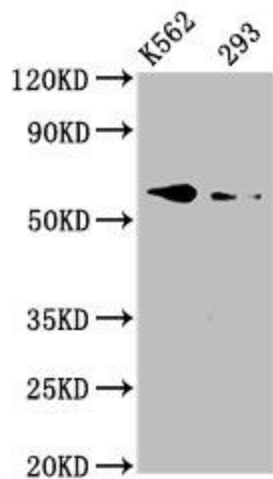
Bone morphogenetic protein receptor type-1A, BMP type-1A receptor, BMPR-1A, Activin receptor-like kinase 3, ALK-3, Serine/threonine-protein kinase receptor R5, SKR5, BMPR1A, ACVRLK3, ALK3

### Application

ELISA, WB, IHC, IF; Recommended dilution: WB:1:500-1:5000, IHC:1:100-1:1000, IF:1:200-1:500



Immunohistochemistry of paraffin-embedded human kidney tissue using BMPR1A Antibody at dilution of 1:100



**Western Blot**

Positive WB detected in: K562 whole cell lysate, 293 whole cell lysate

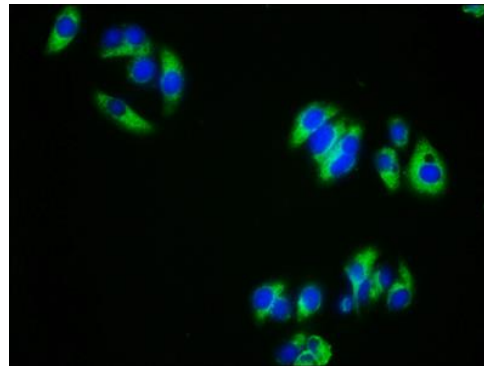
All lanes: BMPR1A antibody at 3.3ug/ml

Secondary

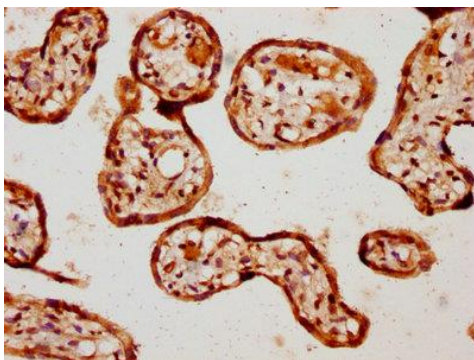
Goat polyclonal to rabbit IgG at 1/50000 dilution

Predicted band size: 61 kDa

Observed band size: 61 kDa



Immunofluorescence staining of HepG2 cells with BMPR1A Antibody at 1:333, counter-stained with DAPI. The cells were fixed in 4% formaldehyde, permeabilized using 0.2% Triton X-100 and blocked in 10% normal Goat Serum. The cells were then incubated with the antibody overnight at 4°C. The secondary antibody was Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L).



IHC image of BMPR1A Antibody diluted at 1:1000 and staining in paraffin-embedded human placenta tissue performed on a Leica Bond™ system. After dewaxing and hydration, antigen retrieval was mediated by high pressure in a citrate buffer (pH 6.0). Section was blocked with 10% normal goat serum 30min at RT. Then primary antibody (1% BSA) was incubated at 4°C overnight. The primary is detected by a biotinylated secondary antibody and visualized using an HRP conjugated SP system.