

BMP7 Polyclonal Antibody

(Catalog # A52444)

Background

Induces cartilage and bone formation. May be the osteoinductive factor responsible for the phenomenon of epithelial osteogenesis. Plays a role in calcium regulation and bone homeostasis.

Description

BMP7 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

P18075

Purification

Protein G purified

Immunogen

Recombinant Human Bone morphogenetic protein 7 protein (293-431AA)

Storage

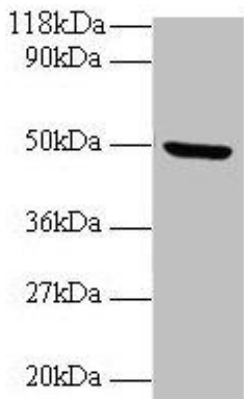
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Osteogenic protein 1, BMP7, OP1

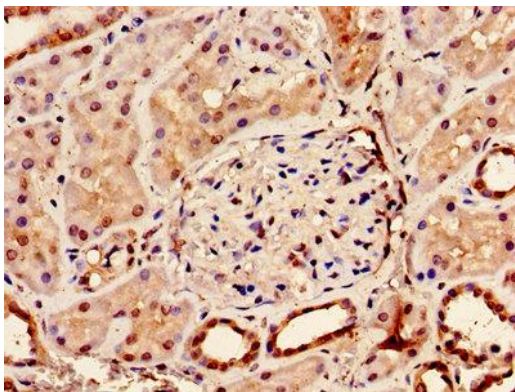
Application

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

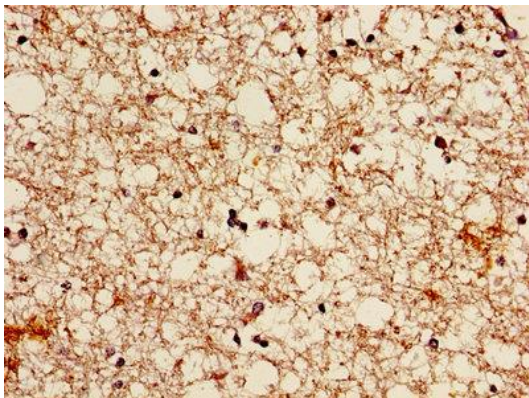


Western blot

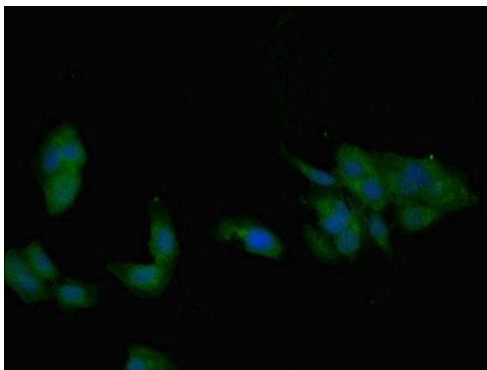
All lanes: BMP7 antibody at 2ug/ml + Mouse kidney tissue
Secondary
Goat polyclonal to rabbit IgG at 1/10000 dilution
Predicted band size: 50 kDa
Observed band size: 50 kDa



Immunohistochemistry of paraffin-embedded human kidney tissue using BMP7 Antibody at dilution of 1:100



Immunohistochemistry of paraffin-embedded human brain tissue using BMP7 Antibody at dilution of 1:100



Immunofluorescent analysis of HeLa cells using BMP7 Antibody at dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)