
BMPER Polyclonal Antibody

(Catalog #A70079)

Background

Inhibitor of bone morphogenetic protein (BMP) function, it may regulate BMP responsiveness of osteoblasts and chondrocytes.

Description

BMPER 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

Q8N8U9

Purification

>95%, Protein G purified

Immunogen

Recombinant Human BMP-binding endothelial regulator protein (271-418AA)

Storage

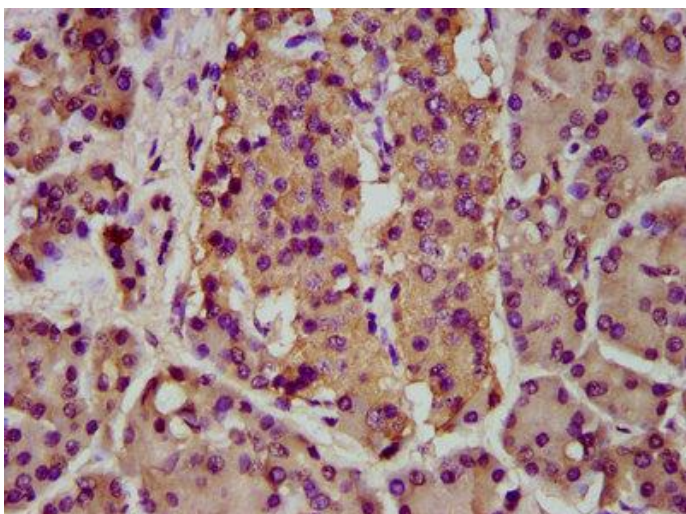
Shipped at 4°C. Upon receipt, store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

BMP-binding endothelial regulator protein, Bone morphogenetic protein-binding endothelial cell precursor-derived regulator, Protein crossveinless-2, hCV2, BMPER, KIAA1965

Application

ELISA, IHC; Recommended dilution: IHC 1:200 - 1:500



IHC image of BMPER Polyclonal Antibody diluted at 1:200 and staining in paraffin-embedded human pancreatic cancer 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.