# Osteopontin polyclonal antibody

Osteopontin (OPN) is an acidic extracellular matrix cell adhesion protein that is relatively abundant not only in bone matrix, plasma, urine, and milk, but is also found in malignant and atherogenic tissues. Phosphorylation, glycosylation and calcium modifications allow intact and fragmented Osteopontin to direct a variety of diverse responses including tissue remodeling, inflammation and cell survival. Plasma Osteopontin has been shown to provide prognostic information in breast, prostate, colon, and lung cancers as well as metastatic carcinomas. The notable presence of Osteopontin in a variety of tumors is strongly correlated to pathological stage, suggesting its critical role in tumor invasiveness, progression and metastasis. In addition, Osteopontin inhibits inducible nitric oxide synthase activity, thereby protecting tumor cells from NO-mediated macrophage cytotoxic attack. Osteopontin is found in atherosclerotic plaques and may drive a number of diabetic vascular pathologies.

This antibody is covered by our Worry-Free Guarantee.

Citations: 2

View Online »

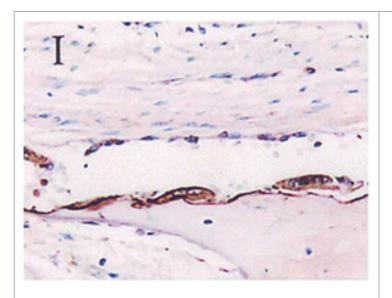
# **Ordering Information**

**Order Online** »

ALX-210-309-R020	20μΙ
ALX-210-309-R100	100μΙ

Manuals, SDS & CofA

View Online »



Immunohistochemistry analysis of fetal calvaria (paraffin section) stained with Osteopontin, pAb (Prod. no. ALX-210-309) at 1:1500.

### **Handling & Storage**

**Handling** Avoid freeze/thaw cycles. After reconstitution, prepare aliquots and store at -20°C.

Long Term Storage +4°C

**Shipping** Blue Ice

## Regulatory Status RUO - Research Use Only

### **Product Details**

Alternative Name Bone sialoprotein 1, Secreted phosphoprotein 1, OPN,

SPP1

Application ELISA, IHC (PS), WB

Crossreactivity Does not cross-react with human osteonectin or bone

sialoprotein.

**Formulation** Lyophilized. From neat/unpurified rabbit serum.

Gene/Protein Identifier 6696 (Entrez GenelD)

**Host** Rabbit

Immunogen Synthetic peptide corresponding to aa 59-74

(P59SKSNESHDHMDDMDD74) of mature human

osteopontin.

Recommendation Dilutions/Conditions ELISA (1:1,000)Immunohistochemistry (paraffin sections,

1:100-1:1500)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be

determined individually for each application.

Reconstitute with 20µl or 100µl distilled water.

Species Reactivity Human, Mouse

UniProt ID P10451

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

•

Last modified: May 29, 2024



info-

eu@enzolifesciences.com