

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.

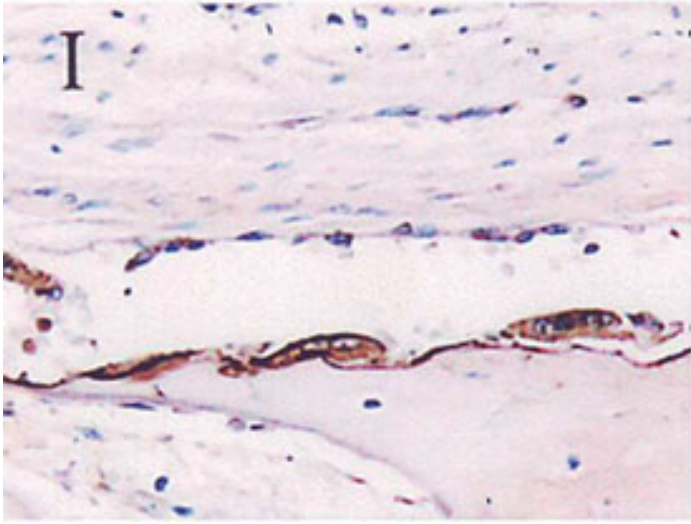
This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 2

[View Online »](#)

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 GeneID)
Host	Rabbit
Immunogen	Synthetic peptide corresponding to aa 59-74 (P59SKSNESHDMDDMDD74) 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.
Reconstitution	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



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)