

# Osteopontin monoclonal antibody (87-B)

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](#).

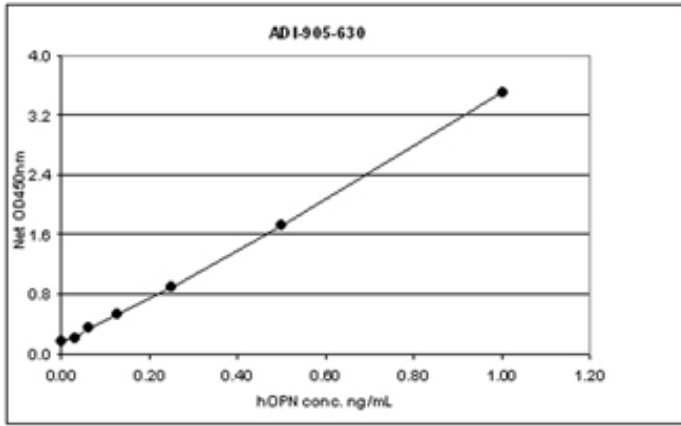
## Ordering Information

[Order Online »](#)

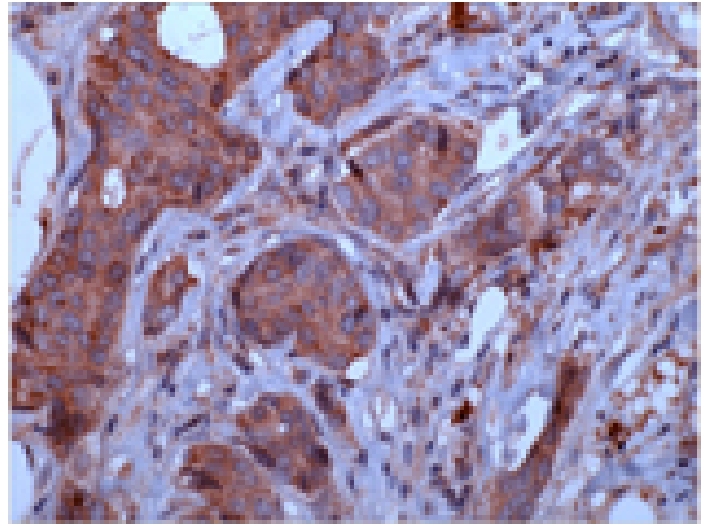
|                 |       |
|-----------------|-------|
| ADI-905-630-100 | 100µg |
|-----------------|-------|

## Manuals, SDS & CofA

[View Online »](#)



Sandwich ELISA Analysis of Osteopontin, mAb (87-B)  
(Prod. No. ADI-905-630)



Immunohistochemistry analysis of formalin-fixed,  
paraffin-embedded breast cancer tissue stained with  
2µg/ml osteopontin, mAb.

## Handling & Storage

**Handling** Avoid freeze/thaw cycles.

**Long Term Storage** -20°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** OPN, SPP1, Secreted phosphoprotein 1, Bone sialoprotein 1

**Application** ELISA, IHC (FS), IHC (PS), WB

**Clone** 87-B

**Crossreactivity** Does not cross-react with mouse or bovine OPN.

**Formulation** Liquid. In PBS, pH 7.2, containing 0.09% sodium azide.

**Host** Mouse

**Immunogen** Recombinant GST tagged human osteopontin.

**Isotype** IgG2a

**Purity Detail** Protein G affinity purified.

**Recommendation Dilutions/Conditions** Immunohistochemistry: staining of formalin-fixed tissues requires boiling tissue sections in 10mM citrate, pH 6.0, for 10min followed by cooling at RT for 20min. Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.

**Source** Purified from mouse ascites.

**Species Reactivity** Human

**Specificity** Recognizes full length OPN and C-terminal portion of thrombin-cleaved OPN.

|                                       |   |
|---------------------------------------|---|
| <b>Technical Info / Product Notes</b> | For immunohistochemistry, staining of formalin-fixed tissues requires boiling tissue sections in 10mM citrate, pH6.0 for 10min followed by cooling at RT for 20min. Recognizes full length human OPN and C-terminal portion of thrombin-cleaved human OPN. Does not react with mouse or bovine OPN. |
| <b>UniProt ID</b>                     | P10451  |
| <b>Worry-free Guarantee</b>           | This antibody is covered by our <a href="#">Worry-Free Guarantee</a> .  |



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)