

# Osteopontin monoclonal antibody (53)

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](#).

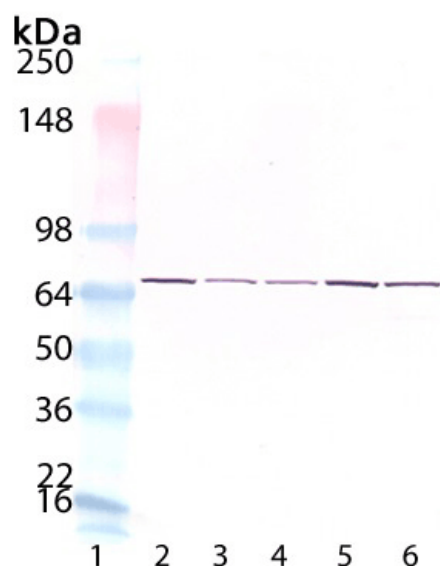
## Ordering Information

[Order Online »](#)

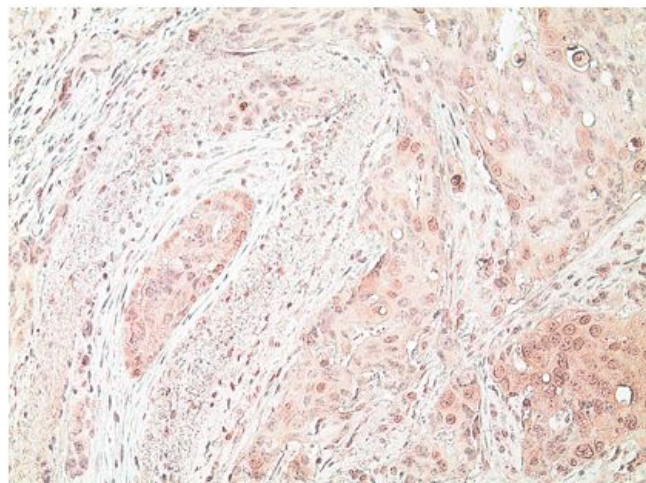
ENZ-ABS753-0100	100µg
-----------------	-------

## Manuals, SDS & CofA

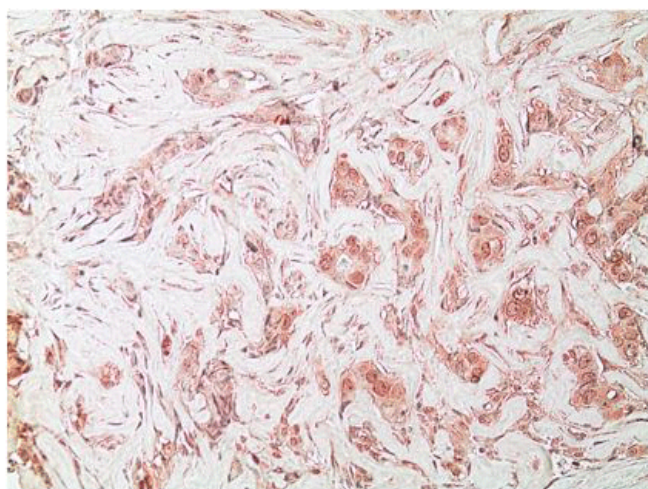
[View Online »](#)



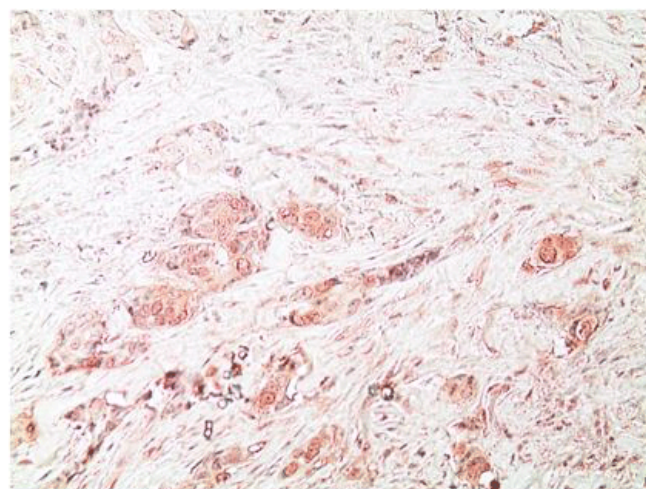
Western Blot analysis of ENZ-ABS753: Lane 1: MW Marker, Lane 2: HeLa lysate, Lane 3: MCF-7 lysate, Lane 4: A431 lysate, Lane 5: HEK-293 lysate, Lane 6: Caki-2 lysate.



Immunohistochemistry analysis of formalin-fixed, paraffin-embedded human breast cancer tissue stained with Osteopontin monoclonal antibody (53), (Prod. No. ENZ-ABS753) at 10 $\mu$ g/ml.



Immunohistochemistry analysis of formalin-fixed, paraffin-embedded human breast cancer tissue stained with Osteopontin monoclonal antibody (53), (Prod. No. ENZ-ABS753) at 10 $\mu$ g/ml.



Immunohistochemistry analysis of formalin-fixed, paraffin-embedded human breast cancer tissue stained with Osteopontin monoclonal antibody (53), (Prod. No. ENZ-ABS753) at 10 $\mu$ g/ml.

<b>Handling</b>	Avoid freeze/thaw cycles.
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Blue Ice

## Product Details

Alternative Name	OPN, SPP1, Secreted phosphoprotein 1, Bone sialoprotein 1
Application	ELISA, IHC (FS), IHC (PS), WB
Clone	53
Formulation	Liquid. In PBS, pH 7.2,containing 0.1% sodium azide.
GenBank ID	U20758
Host	Mouse
Immunogen	Recombinant human Osteopontin.
Isotype	IgG2a
Purity Detail	Protein G affinity purified.
Recommendation Dilutions/Conditions	Western Blot (1:1,000, colorimetric)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Source	Purified from hybridoma tissue culture supernatant.
Species Reactivity	Human
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)