

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](#).

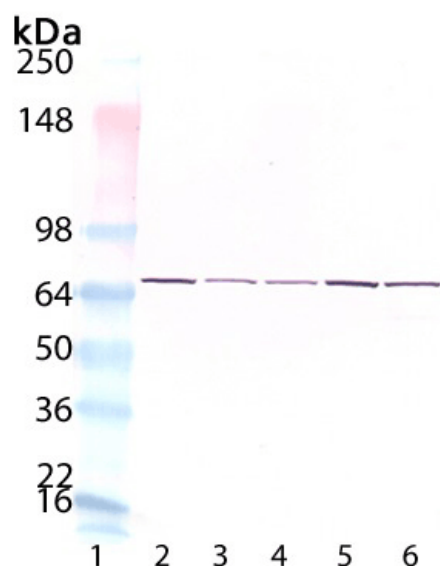
Ordering Information

[Order Online »](#)

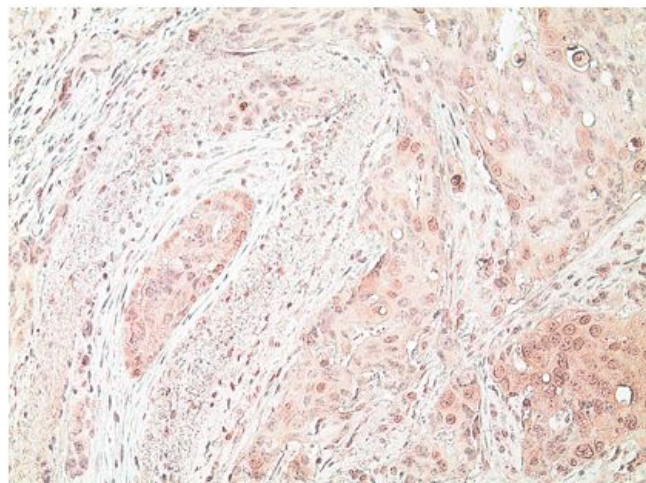
ENZ-ABS753-0100	100µg
-----------------	-------

Manuals, SDS & CofA

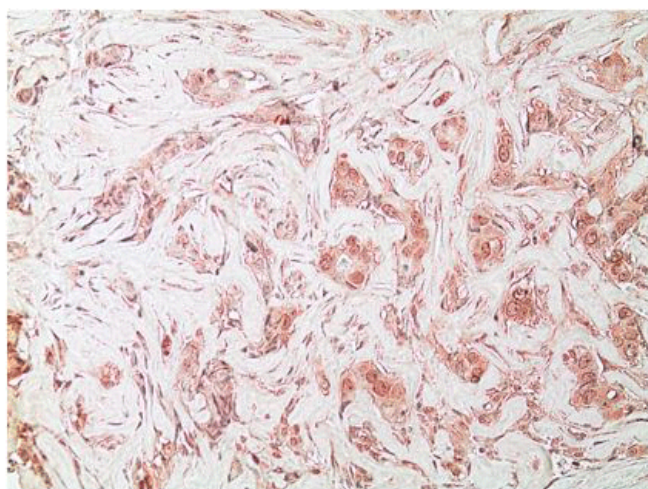
[View Online »](#)



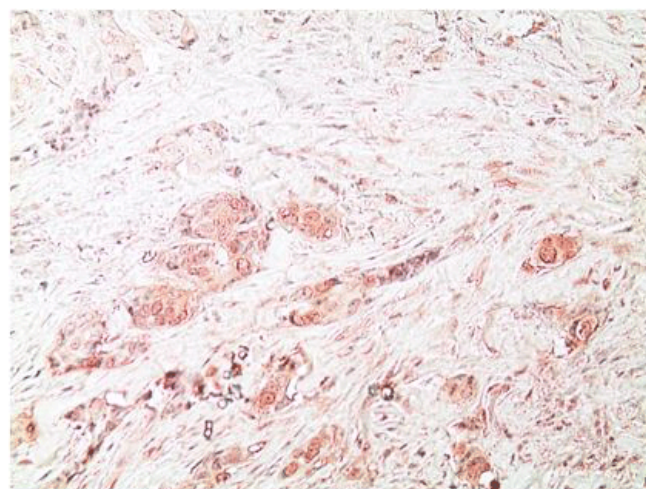
Western Blot analysis of ENZ-ABS753: Lane 1: MW Marker, Lane 2: HeLa lysate, Lane 3: MCF-7 lysate, Lane 4: A431 lysate, Lane 5: HEK-293 lysate, Lane 6: Caki-2 lysate.



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



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



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

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 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 Western Blot (1:1,000, colorimetric) Suggested dilutions/conditions may not be available
Dilutions/Conditions 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



ENZO LIFE SCIENCES,
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 1 402 44 555
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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

Worry Free Guarantee This antibody is covered by our **Worry Free Guarantee**