

Osteopontin recombinant monoclonal antibody (001)

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

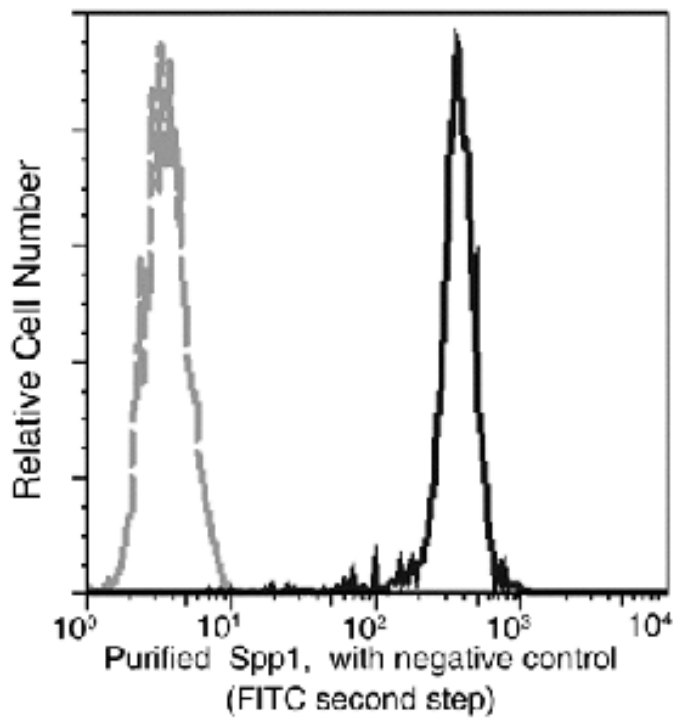
Ordering Information

[Order Online »](#)

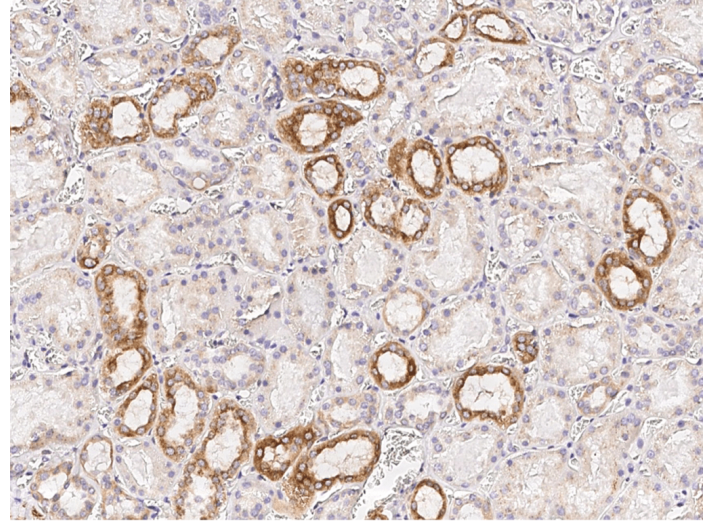
ENZ-ABS944-0100	100µl
-----------------	-------

Manuals, SDS & CofA

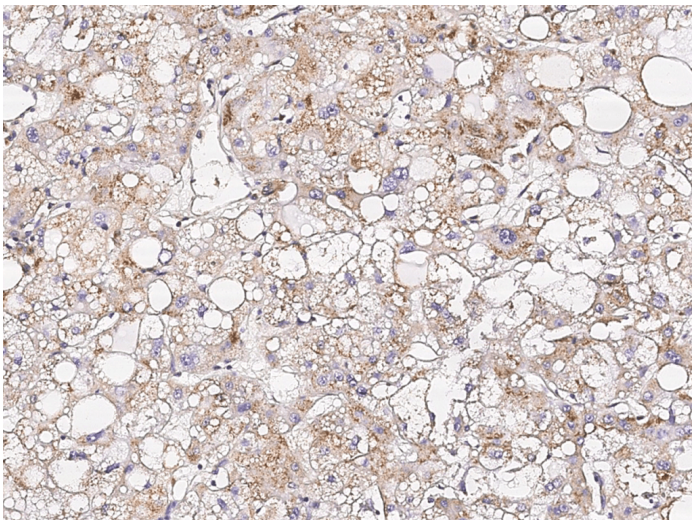
[View Online »](#)



Flow cytometric analysis of purified anti-human Spp1 on U937 cells. U937 cells were treated and stained with Purified Rabbit anti-Spp1 (Bold line histogram), To demonstrate specificity of staining the binding of anti-human Spp1 was blocked by the preincubation of the purified antibody with molar excess of recombinant human Spp1 (5 μ g) for 1 hour (dashed line histogram), then stained with a FITC-conjugated second step antibody.



Immunohistochemical staining of human SPP1 in human kidney with rabbit monoclonal antibody (1:2000, formalin-fixed paraffin embedded sections).



Immunohistochemical staining of human SPP1 in human hepatoma with rabbit monoclonal antibody (1:2000, formalin-fixed paraffin embedded sections).

Handling & Storage

Handling Protected from light. 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, Flow Cytometry, IHC (PS)
Clone	001
Formulation	Liquid. 0.2 µm filtered solution in PBS. Preservative free.
Host	Rabbit
Immunogen	Recombinant human Osteopontin / SPP1 / ETA-1 protein.
Isotype	IgG
Purity Detail	Protein A purified.
Recommendation Dilutions/Conditions	ELISA (1:5'000-1:10'000)Immunohistochemistry (paraffin sections) (1:1'000-1:4000)Flow Cytometry (1:25-1:100)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee



ENZO LIFE SCIENCES,
INC.
Phone: 800 612 0430
[info-](mailto:info-usa@enzolifesciences.com)
usa@enzolifesciences.com

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

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

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

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

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