

TIMP-2 monoclonal antibody (3A4)

TIMP-2 is an inhibitor of matrix metalloproteinases (MMPs) that regulates extracellular matrix remodeling. It is involved in tissue homeostasis, angiogenesis, and cancer progression.

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

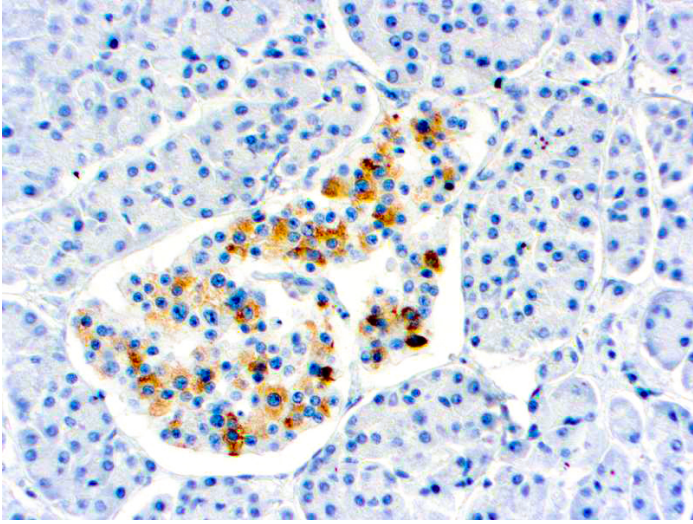
Ordering Information

[Order Online »](#)

ENZ-ABS1089-0500	500µl
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Formalin fixed paraffin embedded human pancreas
stained with TIMP-2 antibody

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Tissue inhibitor of metalloproteinases 2

Application IHC (PS)

Application Notes Pretreatment – EDTA Buffer pH 8.0. Incubation & Temperature 30 min @ RT

Clone 3A4

Formulation Tissue culture supernatant. Contains sodium azide as a preservative.

Host Mouse

Immunogen BALB/C mice were injected with a synthetic peptide from the N-terminal region of human TIMP-2.

Isotype IgG2ak

Positive Control Breast, Colon, Endometrial, HeLa cells, Normal Placenta, Prostate or Ovarian Carcinoma

Recommendation Immunohistochemistry (paraffin sections) (1:50-1:200)

Dilutions/Conditions 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: 800.942.0430
[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-be@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)