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

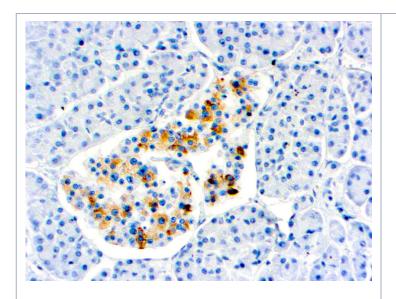
Order Online »

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

ImmunogenBALB/C mice were injected with a synthetic peptide from

the N-terminal region of human TIMP-2.

IgG2ak

Positive Control Breast, Colon, Endometrial, HeLa cells, Normal Placenta,

Prostate or Ovarian Carcinoma

Recommendation Dilutions/Conditions Immunohistochemistry (paraffin sections) (1:50-1:200)

Suggested dilutions/conditions may not be available for all

applications.

Optimal conditions must be determined individually for

each application.

Species Reactivity Human

Last modified: June 6, 2025



info-

eu@enzolifesciences.com