

TIMP-1 (human neutrophils)

Specific inhibitor of matrix metalloproteinases (MMPs). It has no activity on mammalian membrane metalloproteinases, procollagen peptidase or on bacterial metalloproteinases such as thermolysin.

Citations: 6

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-200-426-C005	5µg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. TIMP-1 is very stable if stored at -80°C. It can be kept at -20°C for several months or +4°C for weeks without significant loss of activity.
Handling	Avoid freeze/thaw cycles.
Short Term Storage	-20°C
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Tissue inhibitor of metalloproteinase 1
Formulation	Liquid. In 50mM TRIS-HCl, pH 7.0, containing 200mM sodium chloride, 5mM CaCl ₂ , 1μM ZnCl ₂ , 0.05% BRIJ-35 and 0.05% sodium azide.
MW	~28kDa.
Purity	≥92% (SDS-PAGE, Western blot)
Source	Isolated from stimulated human neutrophils. The secreted protein consists of 184 amino acids, six disulfide bonds and two glycosylation sites containing N-linked oligosaccharides.
Technical Info / Product Notes	Note: It is recommended to perform a preincubation of ~20 min. at 37°C to allow formation of the enzyme-inhibitor complex.
UniProt ID	P01033



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)