

# 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 <sub>2</sub> , 1μM ZnCl <sub>2</sub> , 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	<b>Note:</b> It is recommended to perform a preincubation of ~20 min. at 37°C to allow formation of the enzyme-inhibitor complex.
UniProt ID	P01033

Last modified: May 29, 2024



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-  
fr@enzolifesciences.com](mailto:info-fr@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)