

# TIMP-1 (human) ELISA kit

**Highly sensitive TIMP-1 ELISA kit enabling detection of TIMP-1 in serum, plasma, saliva, urine and tissue culture media in just 2.5 hours.**

The TIMP-1 ELISA kit is a colorimetric immunometric immunoassay kit with results in 2.5 hours.

Tissue inhibitor of metalloproteinases-1, TIMP-1, is one of the four members of the TIMP family which inhibit the activity of the matrix metalloproteinases (MMPs), a large group of peptidases involved in degradation of the extracellular matrix. TIMP-1 displays multiple biological functions. TIMP-1 is capable of inhibiting the activities of most MMPs except membrane-type MMP subfamily. Associated with its inhibitory activity on MMPs, TIMP-1 is able to promote cell proliferation and survival in a wide range of cell types. Recent findings have separated the MMP inhibitory activity of TIMP-1 from its growth promoting effect. TIMP-1 exhibits growth factor-like activity and acts as a cell survival factor which may be mediated through ligand-receptor interactions.

The balance of TIMPs/MMPs plays a crucial role in extracellular matrix homeostasis and remodeling, the imbalance of TIMPs/MMPs is involved in various kinds of diseases in multiple organs. Breaking the balance between MMPs and TIMPs occurs in various pathologic processes and involves all the members of these two families. Elevation of TIMP-1 has been reported in diseases such as systemic sclerosis, amyotrophic lateral sclerosis (ALS), renal disease and invasive tumors. Down regulation of TIMP-1 was found to be associated with ischemic heart failure and chronic inflammation.

## Ordering Information

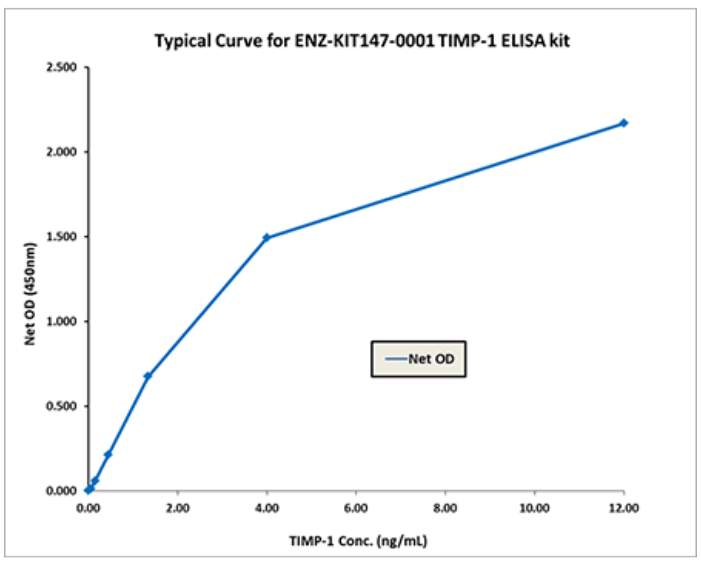
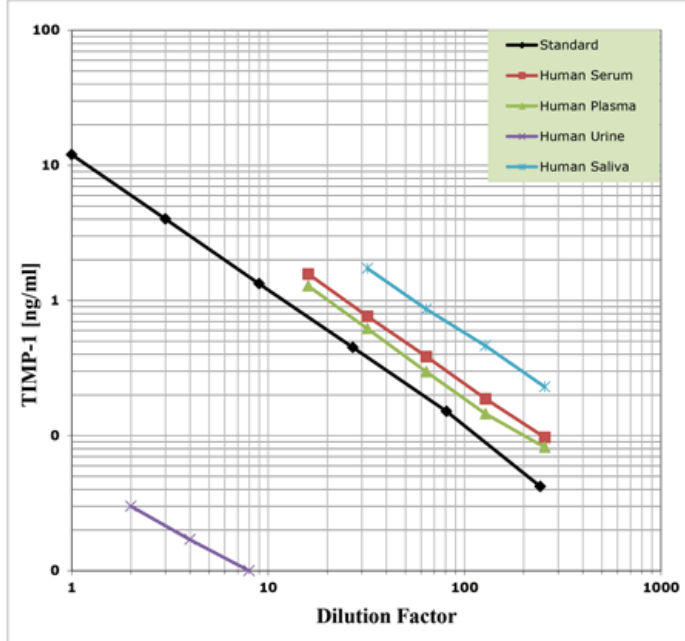
[Order Online »](#)

ENZ-KIT147-0001	96 wells
-----------------	----------

## Manuals, SDS & CofA

[View Online »](#)

- Highly sensitive measurement of TIMP-1, detecting as little as 30 pg/ml
- Negligible cross-reactivity with similar proteins (TIMP-2, TIMP-3, TIMP-4, MMP-1, MMP-2, MMP-3, MMP-7, MMP-8, MMP-9, MMP-13)
- High throughput format with results in 2.5 hours for up to 39 samples in duplicate
- Fully quantitative results that surpass semi-quantitative Western blot analysis



## Handling & Storage

Long Term Storage      -20°C

Shipping      Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	Tissue inhibitor of metalloproteinases-1
Application	Colorimetric detection, ELISA
Application Notes	For the quantitative determination of TIMP-1 in human serum, plasma, saliva, urine and tissue culture media.
Assay Time	2.5 hours
Compatibility	This product is compatible with the <a href="#">Absorbance 96 Plate Reader</a> .
Contents	Microtiter plate, Assay Buffer 13, Standard, Wash Buffer Concentrate, Antibody, Antibody Diluent, Conjugate, Substrate, and Stop Solution
Crossreactivity	Negligible cross-reactivity with similar proteins (TIMP-2, TIMP-3, TIMP-4, MMP-1, MMP-2, MMP-3, MMP-7, MMP-8, MMP-9, MMP-13).
Sensitivity	≤ 30 pg/mL (0.049 – 12 ng/ml)
Species Reactivity	Human
UniProt ID	P01033
Wavelength	450 nm

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)