

# MMP-14 (catalytic domain) (human), (recombinant)

Citations: 7

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-SE259-0010	10µg
----------------	------

Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

<b>Use/Stability</b>	The enzyme is stable on ice for at least several hours. However, it is recommended that thawing and dilution of the enzyme be done within as short a time as possible before start of the assay. NOTE: When stored at -70°C, this enzyme is stable at the concentration supplied, in its current storage buffer. Procedures such as dilution of the enzyme followed by refreezing could lead to loss of activity.
<b>Handling</b>	After opening, prepare aliquots and store at -70°C. Avoid freeze/thaw cycles.
<b>Long Term Storage</b>	-80°C
<b>Shipping</b>	Dry Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

<b>Activity</b>	Preincubation of MMP-14 catalytic domain at 13.6nM with the broad-spectrum inhibitor GM6001 (Prod. No. BML-EI300) at 20nM for 1 hour completely inhibits enzymatic activity.
<b>Alternative Name</b>	Matrix metalloproteinase 14, Membrane type matrix metalloproteinase 1, MT-1 MMP
<b>Application Notes</b>	Useful tool to study enzyme kinetics, cleave target substrates, and screen for inhibitors.
<b>Formulation</b>	Liquid. In 50mM TRIS, pH 7.5, containing 5mM calcium chloride, 300mM sodium chloride, 20µM zinc chloride, 0.5% Brij-35, and 30% glycerol.
<b>MW</b>	22.5 kDa
<b>Purity Detail</b>	Partially purified by single-step affinity chromatography and gel filtration.
<b>Source</b>	Produced in <i>E.coli</i> . Active Matrix Metalloproteinase-14 (MMP-14, Membrane-Type Matrix Metalloproteinase 1, MT1-MMP) catalytic domain from human cDNA. The enzyme consists of the catalytic domain of human MMP-14 (Tyr <sup>112</sup> -Arg <sup>298</sup> , NM_004995) with a C-terminal purification tag. This represents a naturally-occurring active form of MMP-14 which lacks the C-terminal hemopexin domain. MMPs lacking this domain cannot cleave native collagens; however, activity toward other targets such as gelatin, casein, or peptide substrates is unaffected.



ENZO LIFE SCIENCES  
INC.  
Phone: 800.942.0430  
info-  
[usa@enzolifesciences.com](mailto:usa@enzolifesciences.com)

Specific Activity  $\geq 1000$  pmol/min/µg at 37°C using the colorimetric thiopeptide Ac-Pro-Leu-Gly-S-Leu-Leu-Gly-OMe (100 µM; Prod. No. BML-P125) as substrate.

ENZO LIFE SCIENCES  
EUROPEAN DIVISION  
& Luxembourg  
Phone: +41 61 926 8989  
info-  
[eu@enzolifesciences.com](mailto:eu@enzolifesciences.com)

ENZO LIFE SCIENCES  
FRANCE  
Phone: +33 472 440 655  
info-  
[fr@enzolifesciences.com](mailto:fr@enzolifesciences.com)

ENZO LIFE SCIENCES  
GERMANY  
Phone: +49 7621 5500 526  
info-  
[de@enzolifesciences.com](mailto:de@enzolifesciences.com)

ENZO LIFE SCIENCES  
UK  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
info-  
[uk@enzolifesciences.com](mailto:uk@enzolifesciences.com)

UniProt ID

P50281

For local distributors and detailed product information visit us online: [enzo.com](http://enzo.com) All product names, logos, and brands are the property of their respective owners.