

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

Citations: 8

[View Online »](#)

Ordering Information

[Order Online »](#)

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

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

| | |
|-------------------|---|
| Use/Stability | 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. |
| Handling | After opening, prepare aliquots and store at -70°C. Avoid freeze/thaw cycles. |
| Long Term Storage | -80°C |
| Shipping | Dry Ice |

Regulatory Status

RUO - Research Use Only

Product Details

| | |
|-------------------|--|
| Activity | 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. |
| Alternative Name | Matrix metalloproteinase 14, Membrane type matrix metalloproteinase 1, MT-1 MMP |
| Application Notes | Useful tool to study enzyme kinetics, cleave target substrates, and screen for inhibitors. |
| Formulation | Liquid. In 50mM TRIS, pH 7.5, containing 5mM calcium chloride, 300mM sodium chloride, 20µM zinc chloride, 0.5% Brij-35, and 30% glycerol. |
| MW | 22.5 kDa |
| Purity Detail | Partially purified by single-step affinity chromatography and gel filtration. |

| | |
|--------------------------|--|
| Source | 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 ¹¹² -Arg ²⁹⁸ , 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. |
| Specific Activity | ≥1000 pmol/min/μg at 37°C using the colorimetric thiopeptolide Ac-Pro-Leu-Gly-S-Leu-Leu-Gly-OEt (100 μM; Prod. No. BML-P125) as substrate. |
| UniProt ID | P50281 |

Last modified: June 16, 2025



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)