

MMP-8 (catalytic domain) (human), (recombinant)

Citations: 5

[View Online »](#)

Ordering Information

[Order Online »](#)

| | |
|----------------|------|
| BML-SE255-0010 | 10µg |
|----------------|------|

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

| | |
|-------------------|--|
| Activity | Preincubation of MMP-8 catalytic domain at 13.6 nM with the broad-spectrum inhibitor GM6001 (Prod #BML-EI300) at 20nM for 1 hour completely inhibits enzymatic activity. |
| Alternative Name | Matrix metalloproteinase 8, Neutrophil collagenase, Collagenase-2 |
| Application Notes | Useful tool to study enzyme kinetics, cleave target substrates, and screen for inhibitors. |
| Formulation | Liquid. In 50mM TRIS, 5mM CaCl ₂ , 300mM NaCl, 20μM ZnCl ₂ , 0.5% Brij-35, and 30% glycerol. |
| MW | 20.3 kDa |
| Purity | ≥90% (SDS-PAGE) |
| Purity Detail | Purified by multi-step chromatography. |
| Source | Produced in <i>E. coli</i> . Active Matrix Metalloproteinase-8 (MMP-8, neutrophil collagenase, collagenase-2) catalytic domain from human cDNA. The enzyme consists of the catalytic domain of human MMP-8 (Phe ⁹⁹ -Gln ²⁶⁹ , NM_002424) with a C-terminal purification tag. This represents a naturally-occurring active form of MMP-8 which lacks the C-terminal hemopexin domain ¹ . MMPs lacking this domain cannot cleave native collagens; however, activity toward other targets such as gelatin, casein, or peptide substrates is unaffected. |
| Specific Activity | ≥2000 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 | P22894 |



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)