

# 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<sub>2</sub>, 300mM NaCl, 20µM ZnCl<sub>2</sub>, 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 *E. coli*. 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<sup>99</sup>-Gln<sup>269</sup>, NM\_002424) with a C-terminal purification tag. This represents a naturally-occurring active form of MMP-8 which lacks the C-terminal hemopexin domain1. 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)