

# MMP-11 (catalytic domain) (human), (recombinant)

MMP-11 may be a link between obesity and cancer.

Citations: 1

[View Online »](#)

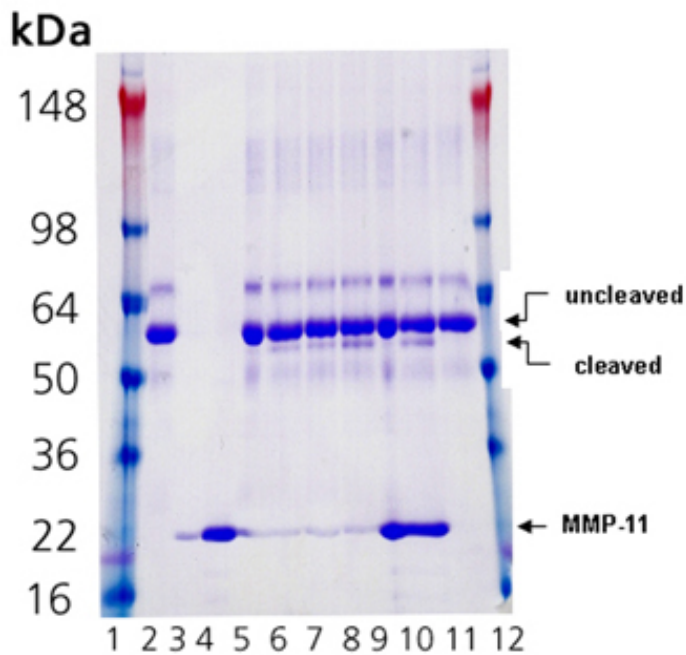
## Ordering Information

[Order Online »](#)

BML-SE282-0010	10µg
----------------	------

Manuals, SDS & CofA

[View Online »](#)



Coomassie-stained SDS-PAGE showing a timecourse of  $\alpha$ 1-antitrypsin (3 $\mu$ g) cleavage after incubation at 37°C with or without MMP-11 catalytic domain. Lanes are as follows: 1. Marker; 2. 1.5  $\mu$ g  $\alpha$ 1-antitrypsin (0 hours); 3. 50 ng MMP-11, (0 hours); 4. 1  $\mu$ g MMP-11, (0 hours); 5. 1.5  $\mu$ g  $\alpha$ 1-antitrypsin incubated with 50 ng MMP-11 (0 hours); 6. 1.5  $\mu$ g  $\alpha$ 1-antitrypsin incubated with 50 ng MMP-11 (3 hours); 7. 1.5  $\mu$ g  $\alpha$ 1-antitrypsin incubated with 50 ng MMP-11 (8 hours); 8. 1.5  $\mu$ g  $\alpha$ 1-antitrypsin incubated with 50 ng MMP-11 (24 hours); 9. 1.5  $\mu$ g  $\alpha$ 1-antitrypsin incubated with 1  $\mu$ g MMP-11 (0 hours); 10. 1.5  $\mu$ g  $\alpha$ 1-antitrypsin incubated with 1  $\mu$ g MMP-11 (24 hours); 11. 1.5  $\mu$ g  $\alpha$ 1-antitrypsin (24 hours); 12. Marker.

## Handling & Storage

Long Term Storage      -80°C

Shipping      Dry Ice

**Regulatory Status**    RUO - Research Use Only

## Product Details

**Alternative Name**      Matrix metalloproteinase 11, Stromelysin-3

**Application Notes**      Useful 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**      19.3 kDa

**Purity**      ≥95% (SDS-PAGE)

**Purity Detail**      Purified by multi-step chromatography.

**Source**      Produced in *E. coli*. Active Matrix Metalloproteinase-11 (MMP-11, Stromelysin-3) catalytic domain from human cDNA. The enzyme consists of the catalytic domain of human MMP-11 (Phe<sup>98</sup>-Ser<sup>266</sup>, NM\_005940) with a C-terminal purification tag. MMPs lacking this domain cannot cleave native collagens; however, activity toward other targets such as gelatin, casein, or peptide substrates is unaffected. It may be an important link between obesity and cancer.

**Specific Activity**      Due to its unusual substrate preferences [A(A/Q)(N/A)~(L/Y)(T/V/M/R)(R/K), or G(G/A)E~LR5], MMP-11 cleaves MMP peptide substrates such as Prod. No. BML-P125, BML-P126, and BML-P132 extremely slowly (several hours yield very little product). Therefore, the activity of each lot of MMP-11 is verified by digestion of macromolecules.

**UniProt ID**      P24347



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)