

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

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

Citations: 1

[View Online »](#)

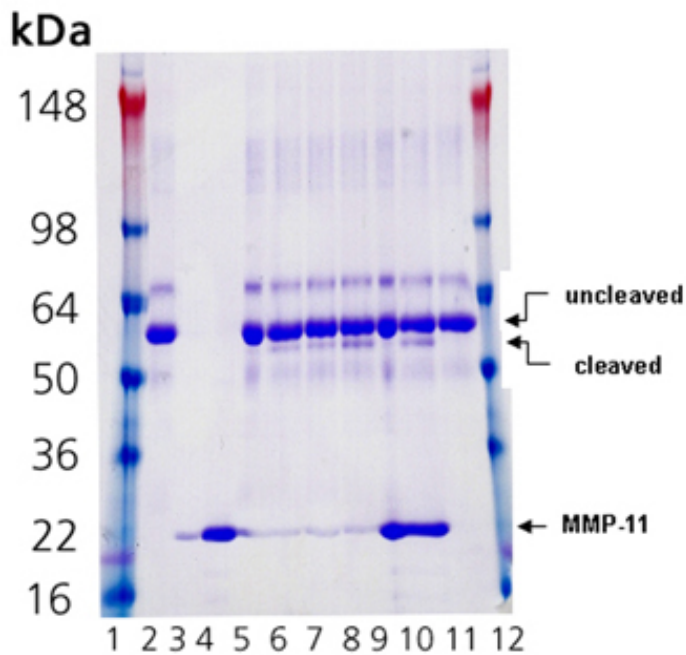
Ordering Information

[Order Online »](#)

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

Manuals, SDS & CofA

[View Online »](#)



Coomassie-stained SDS-PAGE showing a timecourse of α 1-antitrypsin (3 μ g) cleavage after incubation at 37°C with or without MMP-11 catalytic domain. Lanes are as follows: 1. Marker; 2. 1.5 μ g α 1-antitrypsin (0 hours); 3. 50 ng MMP-11, (0 hours); 4. 1 μ g MMP-11, (0 hours); 5. 1.5 μ g α 1-antitrypsin incubated with 50 ng MMP-11 (0 hours); 6. 1.5 μ g α 1-antitrypsin incubated with 50 ng MMP-11 (3 hours); 7. 1.5 μ g α 1-antitrypsin incubated with 50 ng MMP-11 (8 hours); 8. 1.5 μ g α 1-antitrypsin incubated with 50 ng MMP-11 (24 hours); 9. 1.5 μ g α 1-antitrypsin incubated with 1 μ g MMP-11 (0 hours); 10. 1.5 μ g α 1-antitrypsin incubated with 1 μ g MMP-11 (24 hours); 11. 1.5 μ g α 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 CaCl₂, 300mM NaCl, 20μM ZnCl₂, 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 (Phe⁹⁸-Ser²⁶⁶, 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)