

MMP-19 (catalytic domain) (human), (recombinant)

MMP-19 is expressed widely, and degrades aggrecan, fibronectin, type I gelatin, and basement membrane components such as laminin, nidogen, and type IV collagen. It is involved in angiogenesis, cancer, rheumatoid arthritis, multiple sclerosis, and psoriasis.

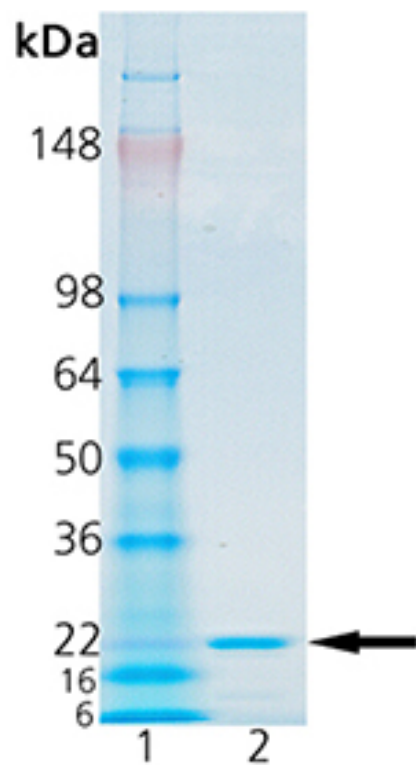
Ordering Information

[Order Online »](#)

BML-SE561-0010	10µg
----------------	------

Manuals, SDS & CofA

[View Online »](#)



SDS-PAGE Analysis: Lane 1: MW Marker, Lane 2: 0.5 μ g of Purified MMP-19 (catalytic domain) (human) (rec.) Prod. No. BML-SE561.

Handling & Storage

Use/Stability	When stored under the recommended conditions, this enzyme is stable at the concentration supplied, in its current storage buffer. Procedures such as dilution followed by refreezing could lead to loss of activity.
Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -80°C.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Activity	Preincubation of MMP-19 catalytic domain at 1.9µM with the broad-spectrum inhibitor GM6001 (Prod No. BML-EI300) at 100nM for 30 minutes only weakly inhibits enzymatic activity (27%).
Alternative Name	Matrix metalloproteinase 19, RASI
Application Notes	Useful tool to study enzyme kinetics, cleave target substrates, and screen for inhibitors.
Formulation	Liquid. In 25mM NaBorate, pH 9.3, containing 5mM CaCl ₂ , 20µM ZnCl ₂ , 0.05% Brij-35 and 20% glycerol.
MW	19.2 kDa
Purity	≥95% (SDS-PAGE)
Purity Detail	Purified by multi-step chromatography.
Sequence	Recombinant MMP-19 corresponding to aa Leu ⁹⁹ -Ser ²⁵⁹ (NM_002429), which comprises the catalytic domain of human MMP-19, with a C-terminal purification tag.
Source	Produced in <i>E.coli</i> . Active recombinant MMP-19.
Specific Activity	≥ 0.40U/µg. One U=100 pmol/min at 37°C using the colorimetric thiopeptolide Ac-Pro-Leu-Gly-S-Leu-Leu-Gly-OEt (100µM; Prod. No. BML-P125) as substrate.



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

Q99542

European Sales Office
ENZO LIFE SCIENCES
(EU) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com