

MMP-13 (catalytic domain) (human), (recombinant)

Citations: 8

[View Online »](#)

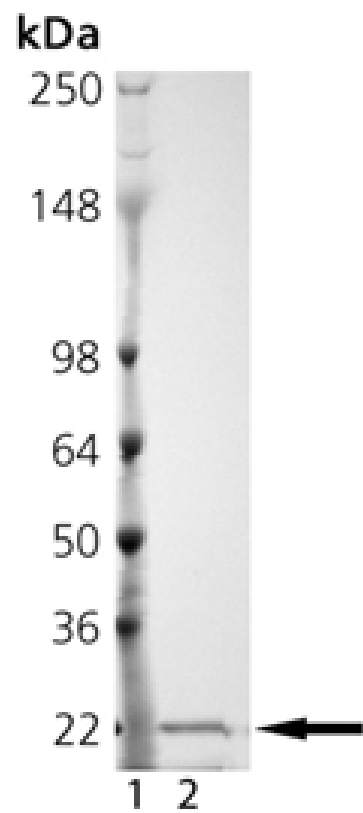
Ordering Information

[Order Online »](#)

BML-SE246-0010	10µg
----------------	------

Manuals, SDS & CofA

[View Online »](#)



SDS-PAGE Analysis: Lane1: MW Marker; Lane 2: 2 μ g of purified MMP-13 (catalytic domain) (human), (recombinant).

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Activity	Preincubation of MMP-13 catalytic domain at 13nM with the broad-spectrum inhibitor GM6001 (Prod. No. BML-EI300) at 20nM for 1 hour completely inhibits enzymatic activity.
Alternative Name	Matrix metalloproteinase 13, Collagenase-3
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	20.4 kDa
Purity	≥95% (SDS-PAGE)
Purity Detail	Purified by multi-step chromatography.
Source	Produced in <i>E. coli</i> . Active Matrix Metalloproteinase-13 (MMP-13, collagenase-3) catalytic domain from human cDNA. The enzyme consists of the catalytic domain of human MMP-13 (Tyr ¹⁰⁴ -Asn ²⁷⁴ , NM_002427) with a C-terminal purification tag. This represents a naturally-occurring active form of MMP-13 which lacks the C-terminal hemopexin domain. 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	P45452



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)