

MMP-10 (catalytic domain) (human), (recombinant)

Citations: 3

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-SE329-0010	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-10 catalytic domain at 22nM with the broad-spectrum inhibitor GM6001 (Prod. No. BML-EI300) at 100nM for 1 hour inhibits enzymatic activity by 95%.
Alternative Name	Matrix metalloproteinase 10, Stromelysin-2
Application Notes	Useful tool to study of 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	19.4 kDa
Purity Detail	Partially purified by single-step affinity chromatography and gel filtration.
Source	Produced in <i>E. coli</i> . Active Matrix Metalloproteinase-10 (MMP-10, stromelysin-2, transin-2) catalytic domain from human cDNA. The enzyme consists of the catalytic domain of human MMP-10 (Phe ⁹⁹ -Glu ²⁷¹ , NM_2425) with a C-terminal purification tag. This comprises an active form of MMP-10 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	≥200 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	P09238



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)