

MMP-3 (catalytic domain) (human), (recombinant)

Citations: 2

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-SE109-0010	10µg
----------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name

Matrix metalloproteinase 3, Stromelysin-1, Transin-1

Application Notes

Study enzyme kinetics, cleave target substrates, and screen for inhibitors.

Formulation

Liquid. In 50mM TRIS, pH 7.5, 5mM CaCl₂, 300mM NaCl, 20μM ZnCl₂, 0.05% Brij-35.

MW

19.5kDa

Purity

≥95% (SDS-PAGE)

Purity Detail

Purified by multi-step chromatography.

Source

Produced in *E. coli*. Active Matrix Metalloproteinase-3 (MMP-3, stromelysin-1, transin) catalytic domain from human cDNA. The enzyme consists of the catalytic domain of human MMP-3 (Phe¹⁰⁰-Thr²⁷², NM_002422) with a C-terminal purification tag.

Specific Activity

≥350 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

P08254

Last modified: May 29, 2024



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)