

MMP-9 (catalytic domain) (human), (recombinant, E. coli)

Matrix metalloproteinase 9 (MMP-9) belongs to a class of enzymes that belong to the zinc-metalloproteinase family involved in the breakdown of extracellular matrix in normal physiological processes, such as embryonic development, reproduction, angiogenesis, bone development, wound healing, cell migration, learning and memory. Its also associated with numerous pathological processes, including cancer, immunologic and cardiovascular diseases.

Citations: 12

[View Online »](#)

Ordering Information

[Order Online »](#)

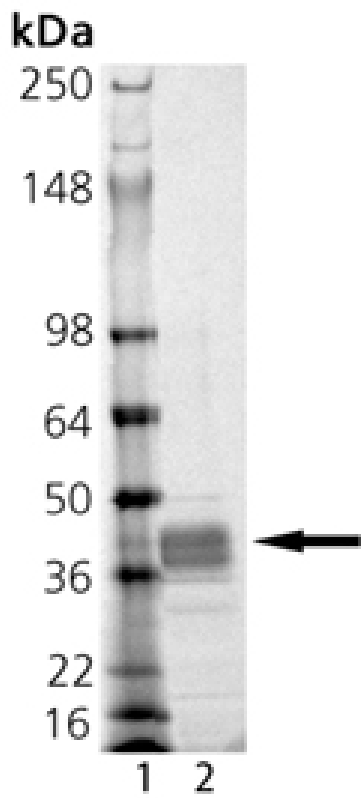
BML-SE360-0010

10µg

Manuals, SDS & CofA

[View Online »](#)

- Naturally-occurring active form of MMP-9
- High purity
- High activity



SDS-PAGE Analysis: Lane 1: MW Marker, Lane 2: 1 μ g MMP-9 (catalytic domain) (human), (recombinant, *E. coli*).

Handling & Storage

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	Yes
Alternative Name	Matrix metalloproteinase 9, Gelatinase B, 92 kDa Type IV collagenase
Application Notes	Useful tool to study enzyme kinetics, cleave target substrates, and screen for inhibitors.
Formulation	Liquid. In 50mM Tris-HCl, pH 7.5, containing 1mM calcium chloride, 300mM sodium chloride, 5µM zinc chloride, 0.1% Brij-35 and 15% glycerol.
MW	39 kDa
Purity	≥95% (SDS-PAGE)
Purity Detail	Purified by multi-step chromatography.
Source	Produced in <i>E. coli</i> . Active recombinant matrix metalloproteinase-9 (MMP-9, gelatinase B, 92 kDa type IV collagenase) cloned from human cDNA. The enzyme consists of residues Phe ¹⁰⁷ -Pro ⁴⁴⁹ (NM_004994), which comprises the catalytic/fibronectin domain of human MMP-9, with a C-terminal purification tag. This represents a naturally-occurring active form of MMP-9 which lacks the C-terminal hemopexin domain. Activity toward its targets, such as gelatin, casein, or peptide substrates, is unaffected.
Specific Activity	≥500 pmol/min/ug at 37°C using the colorimetric thiopeptide Ac-Pro-Leu-Gly-S-Leu-Leu-Gly-OEt (100 µM; Prod. No. BML-P125) as substrate.

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)