

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

BML-SE360

Product Number/Sizes

BML-SE360-0010

10 µg

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

Product Details

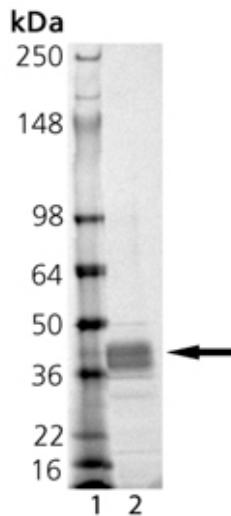
ALTERNATIVE NAME:	Matrix metalloproteinase 9, Gelatinase B, 92 kDa Type IV collagenase
MW:	39 kDa
SOURCE:	Produced in <i>E. coli</i> . Active recombinant matrix metalloproteinase-9 (MMP-9, gelatinase B, 92 kDa type IV collagenase) cloned from human cDNA. The enzyme consists of residues Phe ¹⁰⁷ -Pro ⁴⁴⁹ (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.
UNIPROT ID:	P14780
FORMULATION:	Liquid. In 50mM TRIS, pH 7.5, containing 1mM calcium chloride, 300mM sodium chloride, 5µM zinc chloride, 0.1% Brij-35 and 15% glycerol.
PURITY:	≥95% (SDS-PAGE)
PURITY DETAIL:	Purified by multi-step chromatography.
ACTIVITY:	Yes
SPECIFIC ACTIVITY:	≥500 pmol/min/ug at 37°C using the colorimetric thiopeptolide Ac-Pro-Leu-Gly-S-Leu-Leu-Gly-OEt (100 µM; Prod. No. BML-P125) as substrate.
APPLICATION NOTES:	Useful tool to study enzyme kinetics, cleave target substrates, and screen for inhibitors.
SHIPPING:	Dry Ice
LONG TERM STORAGE:	-80°C
HANDLING:	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -80°C.
SCIENTIFIC BACKGROUND:	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.
REGULATORY STATUS:	RUO - Research Use Only

GLOBAL HEADQUARTERS

Enzo Life Sciences, Inc.
10 Executive Blvd
Farmingdale, NY 11735
USA
T 1-800-942-0430
T 1-631-694-7070
F 1-610-941-9252
E info-usa@enzolifesciences.com
www.enzolifesciences.com

EUROPE/ASIA

Enzo Life Sciences (ELS) AG
Industriestrasse 17, Postfach
CH-4415 Lausen
Switzerland
T +41/061 926 89 89
F +41/061 926 89 79
E info-ch@enzolifesciences.com
www.enzolifesciences.com



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

Product Literature References

- Mechanical strain modulates extracellular matrix degradation and byproducts in an isoform-specific manner* A. Yeganegi, et al. *Biochim. Biophys. Acta Gen. Subj.* **1867** 130286 (2023)
- Validation of Matrix Metalloproteinase-9 (MMP-9) as a Novel Target for Treatment of Diabetic Foot Ulcers in Humans and Discovery of a Potent and Selective Small-Molecule MMP-9 Inhibitor That Accelerates Healing* T.T. Nguyen, et al. *J. Med. Chem.* **61** 8825 (2021)
- Cell-specific expression of the transcriptional regulator RHAMM provides a timing mechanism that controls appropriate wound re-epithelialization* C. Tolg, et al. *J. Biol. Chem.* **295** 5427 (2020)
- Nucleic acid-induced potentiation of matrix metalloproteinase-9 enzymatic activity* T. Duellman, et al. *Biochem. J.* **475** 1597 (2018)
- Substitutions for arginine at position 780 in triple helical domain of the $\alpha 1(I)$ chain alter folding of the type I procollagen molecule and cause osteogenesis imperfecta* E. Makareeva, et al. *PLoS One* **13** e0200264 (2018)
- dCas9-mediated transcriptional activation of tissue inhibitor of metalloproteinases* T. Duellman, et al. *Metalloproteinases Med.* **4** 63 (2017)
- Toxicological effects of NCKU-21, a phenanthrene derivative, on cell growth and migration of A549 and CL1-5 human lung adenocarcinoma cells* H.F. Liao, et al. *PLoS One* **12** e28945763 (2017)
- Active Matrix Metalloprotease-9 Is Associated with the Collagen Capsule Surrounding the Madurella mycetomatis Grain in Mycetoma* K. Geneugelijk, et al. *PLoS Negl. Trop. Dis.* **8** e2754 (2014)
- Directed evolution of protease beacons that enable sensitive detection of endogenous MT1-MMP activity in tumor cell lines* A. Jabaiah, et al. *Chem. Biol.* **18** 392 (2011)
- IL-1b Is Overexpressed and Aberrantly Regulated in Corticosteroid Nonresponders with Autoimmune Inner Ear Disease* S. Pathak, et al. *J. Immunol.* **186** 1870 (2011)

Revised 19-Dec-23

GLOBAL HEADQUARTERS

Enzo Life Sciences, Inc.
10 Executive Blvd
Farmingdale, NY 11735
USA
T 1-800-942-0430
T 1-631-694-7070
F 1-610-941-9252
E info-usa@enzolifesciences.com
www.enzolifesciences.com

EUROPE/ASIA

Enzo Life Sciences (ELS) AG
Industriestrasse 17, Postfach
CH-4415 Lausen
Switzerland
T +41/061 926 89 89
T +41/061 926 89 79
E info-ch@enzolifesciences.com
www.enzolifesciences.com