

MMP-20 (catalytic domain) (human), (recombinant)

MMP-20 functions in tooth development and is implicated in cancer and oral pathology.

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-SE540-0010	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-20 catalytic domain at 0.4nM with the broad-spectrum inhibitor GM6001 (Prod. No. BML-EI300) at 0.1nM for 1 hour completely inhibits enzymatic activity.
Alternative Name	Matrix metalloproteinase 20, Enamelysin
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	19.9 kDa
Purity	≥95% (SDS-PAGE)
Purity Detail	Purified by multi-step chromatography.
Source	Produced in <i>E. coli</i> . Active recombinant matrix metalloproteinase-20 (MMP-20, enamelysin) cloned from human cDNA. The enzyme consists of residues Tyr ¹⁰⁸ -Phe ²⁷⁶ (NM_004771), which comprises the catalytic domain of human MMP-20, with a C-terminal purification tag. This represents a naturally-occurring active form of MMP-20 which lacks the C-terminal hemopexin domain.
Specific Activity	≥ 10000pmol/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	O60882



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)