

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 CaCl₂, 300mM NaCl, 20μM ZnCl₂, 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 *E. coli*. Active recombinant matrix metalloproteinase-20 (MMP-20, enamelysin) cloned from human cDNA. The enzyme consists of residues Tyr¹⁰⁸-Phe²⁷⁶ (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-
fr@enzolifesciences.com](mailto:info-fr@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)