

MMP-9 (human neutrophils), dimer

MMP-9 hydrolyzes the extracellular matrix, e.g. collagen types IV, V and IX and gelatin.

Citations: 3

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-200-423-C005	5µg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Avoid freeze/thaw cycles.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Matrix metalloproteinase 9, Gelatinase B, 92kDa Type IV collagenase
Formulation	Liquid. In 50mM Tris-HCl, pH 7.0, containing 200mM sodium chloride, 5mM CaCl ₂ , 1μM ZnCl ₂ , 0.05% BRIJ35, and 0.05% sodium azide.
MW	~220kDa.
Purity	≥95% (SDS-PAGE, Western blot)
Source	Isolated from stimulated human neutrophil granulocytes (buffy coat). Disulfide-bridged MMP-9 homodimer.
Specific Activity	≥1,400mU/mg protein. One unit is defined as the amount of enzyme that hydrolyzes 1μmol Dnp-Pro-Gln-Gly-Ile-Ala-Gly-Gln-D-Arg-OH per min. at 37°C, pH 7.0.
UniProt ID	P14780

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com