

MMP-8 proenzyme (human neutrophils)

MMP-8 is a glycoprotein containing complex N-linked oligosaccharides. It hydrolyzes type I over type II, and III collagens. Activated MMP-8 is inhibited by TIMP-1 (Prod. No. ALX-200-426).

Citations: 5

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-200-421-C005	5µg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for 1 week when stored at +4°C and for several weeks when stored at -20°C.
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 8, Neutrophil collagenase, Collagenase-2
Concentration	100µg/ml
Formulation	Liquid. In 50mM TRIS-HCl, pH 7.0, containing 200mM NaCl, 5mM CaCl ₂ , 1µM ZnCl ₂ , 0.05% BRIJ 35, and 0.05% sodium azide.
MW	~85kDa.
Purity	≥90% (SDS-PAGE)
Source	Isolated from stimulated human neutrophil granulocytes (buffy coat). Requires activation.
Specific Activity	≥60mU/mg protein (Y. Masui, et al.; Biochem. Med. 17 , 215 (1977)). One unit is defined as the amount of enzyme that hydrolyzes 1µmol 2,4-dinitrophenyl-Pro-Gln-Gly-Ile-Ala-Gly-Gln-D-Arg-OH per min. at 37°C, pH 7.0.
Technical Info / Product Notes	Precursor enzyme needs activation using 2mM AMPA (aminophenylmercuric acetate) or 1mM mersalyl acid for 60 min. at 37°C. Alternatively use 0.1mM PCMB (p-chloromercuribenzoate) or 10µg/ml trypsin for 20 min. at 25°C; PCMB is substantially more effective.
UniProt ID	P22894



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