

# Collagenase AF-1 GMP Grade from Clostridium histolyticum

Collagenase AF-1 GMP Grade is derived from C.histolyticum using a production process completely free of any animal-based components.

## Ordering Information

[Order Online »](#)

NB-17457-02	150U
-------------	------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Activity** Collagenase  $\geq 150$  U/Vial or  $\geq 2000$  U/Vial (PZ activity according to Wunsch, 25 °C) Collagenase  $\geq 3000$  U/mg lyophilisate (PZ activity according to Wunsch, 25 °C) Neutral Protease  $\leq 0.050$  U/mg lyophilisate (DMC assay) PZ Unit (acc. to Wunsch): 1 U according to Wunsch catalyzes the hydrolysis of 1  $\mu$ mol 4-phenylazobenzoyloxycarbonyl-L-prolyl-L-leucylglycyl-L-prolyl-D-arginine per minute at 25 °C, pH 7.1. Neutral Protease: 1 DMC Unit catalyzes the cleavage of 1  $\mu$ mol peptide bond from dimethylcasein per minute at 25 °C, pH 7.0, expressed in terms of newly formed terminal amino groups, determined with TNBS.

**CAS** 9001-12-1

**Endotoxin Content** Endotoxin\*: Sterility (Ph. Eur.) Absence of Clostridia TAMC (Total Aerobic Microbial Count)  $\leq 10$ /Vial TYMC (Total Yeasts/ Moulds Count)  $\leq 10$ /Vial\* Endotoxin: Ph. Eur. (1 Endotoxin Unit is equal to 1 International Unit of a WHO approved reference standard endotoxin (RSE)).

**MW** 70 – 120 kDa (collagenases)

**Purity Detail** Manufactured free from animal components & according to international GMP guidelines.

**Source** From *Clostridium histolyticum*. Contains Collagenase class I and class II.

**Technical Info / Product** NB-17457-01 corresponds to supplier's Cat. No. N0003554

**Notes** For the Original Manufacturer's data sheet please [click here](#).



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)