

# Collagenase NB 4 Standard Grade from Clostridium histolyticum

Collagenase NB 4 Standard Grade is a natural enzyme mixture that contains collagenolytic and additional enzymatic activities including clostripain and neutral protease. The balanced ratio of these activities ensures gentle and efficient tissue dissociation.

## Ordering Information

[Order Online »](#)

NB-17454-02	500mg
NB-17454-01	1g
NB-17454-03	5g

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage +4°C

Shipping Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**Activity** Collagenase activity:  $\geq 0.10$  U/mg (PZ activity acc. to Wünsch, 25 °C)PZ Unit (acc. to Wünsch):1 U according to Wünsch 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.

**CAS** 9001-12-1

**MW** 70 – 120 kDa (collagenases)

**Source** From *Clostridium histolyticum*. Contains class I and class II collagenase 1 and a balanced ratio of proteolytic side activities.

**Technical Info / Product Notes** NB-17454-02 corresponds to supplier's Cat. No. S1745402  
NB-17454-01 corresponds to supplier's Cat. No. S1745401  
NB-17454-03 corresponds to supplier's Cat. No. S1745403

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)