

# Collagenase NB 8

## Broad Range from

### Clostridium

### histolyticum

Collagenase NB 8 Broad Range is a chromatographically purified enzyme resulting in a high collagenolytic activity. This product has an increased collagenase activity compared to Collagenase NB 4 Standard Grade, Collagenase NB 4G Proved Grade and Collagenase NB 7D Gentle Grade. The additional enzymatic activities like clostripain, trypsin-like activity and neutral protease, as well as the endotoxin level are reduced. Collagenase NB 8 is suitable for a broad range of different applications in tissue dissociation such as isolation of rodent pancreatic islets of Langerhans, human cardiomyocytes, tumor cells, muscle cells, osteoblasts and porcine hepatocytes.

#### Ordering Information

[Order Online »](#)

NB-17456-01	250mg
NB-17456-02	1g

#### 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 activity:  $\geq 0.90$  U/mg (PZ activity\*\* acc. to Wunsch, 25 °C)PZ Unit (acc. to Wunsch): 1 U according to Wunsch catalyzes the hydrolysis of 1  $\mu\text{mol}$  4-phenylazobenzoyloxycarbonyl-L-prolyl-L-leucylglycyl-L-prolyl-D-arginine per minute at 25 °C, pH 7.1.

**CAS** 9001-12-1

**Endotoxin Content** Endotoxin\*:  $\leq 100$  EU/mg\*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** Chromatographically purified.

**Source** From *Clostridium histolyticum*. Contains class I and class II collagenase as well as reduced levels of clostripain, trypsin-like activities and neutral protease.

**Technical Info / Product Notes** NB-17456-01 corresponds to supplier's Cat. No. S1745601  
NB-17456-02 corresponds to supplier's Cat. No. S1745602

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)