

# Collagenase NB 4G

## Proved Grade from

### Clostridium

### histolyticum

Collagenase NB 4G Proved Grade is an enzyme mixture which contains collagenolytic and additional enzymatic activities including clostripain and neutral protease. The ratio of these activities leads to gentle and efficient tissue dissociation. The product is suitable for isolation of a broad variety of cell types, especially for fibroblasts, human and rodent hepatocytes, frog oocytes and epithelial cells.

#### Ordering Information

[Order Online »](#)

NB-17465-01	500mg
NB-17465-02	1g
NB-17465-03	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:  $\geq 0.18$  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 collagenase1 and a balanced mix of neutral protease, clostripain and trypsin-like activities.

**Technical Info / Product Notes** NB-17465-01 corresponds to supplier's Cat. No. S1746501  
NB-17465-02 corresponds to supplier's Cat. No. S1746502  
NB-17465-03 corresponds to supplier's Cat. No. S1746503

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)