

Collagenase NB 5 Sterile Grade from Clostridium histolyticum

Collagenase NB 5 Sterile Grade is a natural enzyme mixture that contains collagenolytic and additional enzymatic activities including clostripain and neutral protease and is sterile according to European Pharmacopoeia. The balanced ratio of these activities ensures gentle and efficient tissue dissociation. Collagenase NB 5 is applicable to tissue dissociation such as isolation of human and rodent hepatocytes, chondrocytes, neuronal and endothelial cells, fibroblasts, cells from adipose tissue (e.g. ADSCs), stem cells and their passaging, and rodent osteoblasts.

Ordering Information

[Order Online »](#)

NB-17459-03	1g
-------------	----

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: ≥ 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 μ 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 <i>Clostridium histolyticum</i> . Contains class I and class II collagenase 1 and a balanced ratio of proteolytic side activities.
Technical Info / Product Notes	NB-17459-03 corresponds to supplier's Cat. No. N0002778 For the Original Manufacturer's data sheet please click here .



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-de@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-uk@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com