

Collagenase NB 6

GMP Grade from

Clostridium

histolyticum

Collagenases NB 6 GMP Grade from Clostridium histolyticum is a natural mixture of enzymatic activities and consists of collagenase class I and class II, neutral protease and clostripain. It is used for tissue dissociation such as isolation of human ADSCs, stem cells, chondrocytes, fibroblasts, neuronal and endothelial cells.

Ordering Information

[Order Online »](#)

NB-17458-04	100mg
NB-17458-03	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: ≥ 0.100 PZ U/mg lyophilisate (PZ activity acc. to Wunsch, 25 °C) PZ Unit (acc. to Wunsch): 1 U according to Wunsch catalyzes the hydrolysis of 1 μ mol 4-phenylazobenzyloxycarbonyl-L-prolyl-L-leucylglycyl-L-prolyl-D-arginine per minute at 25 °C, pH 7.1
CAS	9001-12-1
Endotoxin Content	Endotoxin*: Sterility (Ph. Eur.) Absence of Clostridia* 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)
Source	From <i>Clostridium histolyticum</i> . Contains Collagenase class I and class II as well as proteolytic activities such as neutral protease and clostripain.
Technical Info / Product Notes	NB-17458-03 corresponds to supplier's Cat. No. N0002779 NB-17458-04 corresponds to supplier's Cat. No. N0002880. 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-
fr@enzolifesciences.com](mailto:info-fr@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)