

Collagenase NB 1 Premium Grade from Clostridium histolyticum

Collagenase NB 1 Premium Grade is a highly purified collagenase grade used for isolation of sensitive cells and guarantees high yields and viability.

Ordering Information

[Order Online »](#)

| | |
|-------------|-------|
| NB-17455-03 | 2000U |
|-------------|-------|

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: ≥ 3.00 U/mg lyophilisate (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

Endotoxin Content Endotoxin*: ≤ 10.0 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 highly purified class I and class II collagenase. Largely free from clostripain, trypsin-like protease and neutral protease.

Source From *Clostridium histolyticum*. Contains class I and class II collagenase.

Technical Info / Product NB-17455-03 corresponds to supplier's Cat. No. S1745503

Notes 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-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

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

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