

Collagenase NB 7D Gentle Grade from *Clostridium* *histolyticum*

Collagenase NB 7D Gentle Grade is suitable for isolation of a broad variety of cells. Moreover, it can be used especially where sensitive conditions are favored, such as during isolation of dendritic and tumor cells from human or animal tissues. To meet these requirements, Collagenase NB 7D Gentle Grade combines the benefits of slightly higher collagenase activity with low tryptic activities ideal to preserve cell-surface proteins.

Ordering Information

[Order Online »](#)

NB-S17474-01	500mg
NB-S17474-02	1g
NB-S17474-03	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: ≥ 0.2 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 μmol 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 collagenases and reduced trypsin-like activities.

Technical Info / Product 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