

Neutral Protease AF GMP Grade from *Clostridium* *histolyticum*

Neutral Protease AF GMP Grade is derived from *C.histolyticum* using a production process completely free of any animal-based components.

Ordering Information

[Order Online »](#)

NB-30306-01	1Vial
-------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Activity Neutral Protease activity: ≥ 0.50 U/mg lyophilisate (DMC assay) Neutral Protease: 1 DMC Unit catalyzes the cleavage of 1 μ mole peptide bond from dimethylcasein per minute at 25 °C, pH 7.0, expressed in terms of newly formed terminal amino groups, determined with TNBS.

CAS 9001-92-7

Endotoxin Content Endotoxin*: Sterility (Ph. Eur.) Absence of Clostridia TAMC (Total Aerobic Microbial Count) ≤ 10 /Vial TYMC (Total Yeasts/ Moulds Count) ≤ 10 /Vial* Endotoxin: Ph. Eur. (1 Endotoxin Unit is equal to 1 International Unit of a WHO approved reference standard endotoxin (RSE)).

MW 34 kDa

Purity Detail Manufactured free from animal components & according to international GMP guidelines.

Quantity Vial contains not less than 100 DMC U of neutral protease.

Source From *Clostridium histolyticum*.

Technical Info / Product NB-30306-01 corresponds to supplier's Cat. No. N0003553

Notes 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)