

Neutral Protease NB from Clostridium histolyticum

Neutral Protease NB is a metalloprotease suitable for tissue dissociation applications. This enzyme cuts diverse peptides before hydrophobic amino acids and thereby creates smaller fragments.

Ordering Information

[Order Online »](#)

NB-30301-11	50U
NB-30301-12	100U

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 (DMC assay) Neutral Protease: 1 DMC Unit catalyzes the cleavage of 1 μ mol 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

MW 34 kDa

Purity Detail Chromatographically purified.

Source From *Clostridium histolyticum*.

Technical Info / Product Notes NB-30301-11 corresponds to supplier's Cat. No. S3030111
NB-30301-12 corresponds to supplier's Cat. No. S3030112

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