

Neutral Protease NB High Active Grade from Clostridium histolyticum

Neutral Protease NB High Active Grade is a chromatographically purified metalloprotease suitable for different tissue dissociation applications. This product contains an increased proteolytic activity, is largely free from collagenolytic activity (PZ U) and has a very low endotoxin level.

Ordering Information

[Order Online »](#)

NB-30302-02	80U
-------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Activity	Neutral Protease activity: ≥ 4.00 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
Endotoxin Content	Endotoxin*: ≤ 50 EU/mg*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	Chromatographically purified.
Source	From <i>Clostridium histolyticum</i> .
Technical Info / Product Notes	NB-30302-02 corresponds to supplier's Cat. No. S3030202 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