

Clostripain NB from Clostridium histolyticum

Clostripain also known as Clostridiopeptidase B or Endoproteinase Arg-C is a proteinase that cleaves peptide bonds on the carboxyl side of arginine. This protein is a heterodimer composed of a heavy and a light chain that are held together via strong noncovalent forces. As a proteolytic enzyme it can be used for isolation of various cell types by tissue dissociation.

Clostripain exists in two different forms, an oxidized, native condition and a reduced activated form that differ in enzymatic activity. The native clostripain can be activated with reducing agents (e.g. DTT, Cysteine) resulting in a higher activity of the enzyme.

Ordering Information

[Order Online »](#)

NB-S20205-01	200U
--------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Activity Clostripain, native state activity: ≥ 200 U/Vial (BAEE)*Clostripain, native state activity: ≥ 5.0 U/mg (BAEE)*Clostripain, activated activity: ≥ 50.0 U/mg (BAEE)**Clostripain, native state (BAEE assay)1 Unit catalyzes the cleavage of 1 μ mol N α -Benzoyl-L-arginine ethyl ester per minute at 25 °C, pH 7.8.**Clostripain, activated (BAEE assay)1 Unit catalyzes the cleavage of 1 μ mol N α -Benzoyl-L-arginine ethyl ester per minute at 25 °C, pH 7.8. The enzyme is activated with 2.5 mM DTT.

CAS 9028-00-6

MW 59 kDa

Purity Detail Chromatographically highly purified.

Source From *Clostridium histolyticum*.

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