

N-Acetyl-eglin C, (recombinant)

Protease inhibitor

Eglin C was originally isolated from the leech *Hirudo medicinalis* .
Recombinant N-acetyl-eglin C is a 70 amino acid peptide with identical biological activity to native eglin C. An effective inhibitor of chymotrypsin and subtilisin as well as of leukocyte elastase and cathepsin G.

Citations: 2

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status

 RUO - Research Use Only

Product Details

Appearance	White to off-white powder.
CAS	96380-69-7
Formula	$C_{377}H_{552}N_{96}O_{107}$
MW	8141.1
Purity	≥97% (HPLC)
Sequence	Ac-Thr-Glu-Phe-Gly-Ser-Glu-Leu-Lys-Ser-Phe-Pro-Glu-Val-Val-Gly-Lys-Thr-Val-Asp-Gln-Ala-Arg-Glu-Tyr-Phe-Thr-Leu-His-Tyr-Pro-Gln-Tyr-Asn-Val-Tyr-Phe-Leu-Pro-Glu-Gly-Ser-Pro-Val-Thr-Leu-Asp-Leu-Arg-Tyr-Asp-Arg-Val-Arg-Val-Phe-Tyr-Asn-Pro-Gly-Thr-Asn-Val-Val-Asn-His-Val-Pro-His-Val-Gly-OH
Solubility	Soluble in water.
Source	Produced in <i>E. coli</i> .

Last modified: May 29, 2024



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