

Adenylate cyclase toxin

Increases intracellular cAMP levels

AC toxin inserts itself into the membrane of target cells and delivers its adenylate cyclase enzymatic domain to the cell interior. Its adenylate cyclase activity is then stimulated by endogenous Ca^{2+} and calmodulin. The resulting increase in cAMP concentration can easily reach supraphysiological levels.

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-CN150A-0050	50µg
-----------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	After thawing, to ensure recovery of vial contents, centrifuge before opening the tube. Aseptic handling is recommended, no preservatives have been added to the product. Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	AC toxin
Application Notes	AC toxin is a unique tool for increasing cAMP levels in mammalian cells.
Formulation	Liquid. In 0.05M Tris-HCl and 8M Urea pH 8.0.
MW	~200 kDa
Purity	≥80% (SDS-PAGE)
Source	Produced in <i>E. coli</i> .
UniProt ID	J7QLC0



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)