# 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<sup>2+</sup> 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 E. coli.

UniProt ID J7QLC0

Last modified: May 29, 2024

