

ϵ -cAMP

Fluorescent analogue of cyclic AMP

Fluorescent cAMP analogue. Excitation 300nm, Emission 415nm.

*BLG-E001-50 (5 x 10 μ mol pack size) is not sold in the U.S. or Canada.
Please [contact us](#) for available options.*

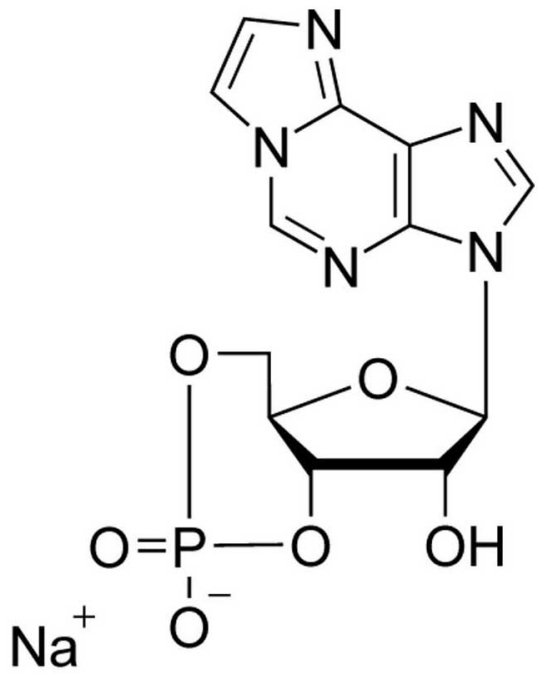
Ordering Information

[Order Online »](#)

| | |
|-------------|----------------|
| BLG-E001-50 | 5x10 μ mol |
|-------------|----------------|

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

| | |
|-------------------|---------------------|
| Handling | Protect from light. |
| Long Term Storage | -20°C |
| Shipping | Ambient Temperature |

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 1,N6-Ethenoadenosine-3',5'-cyclic monophosphate . sodium salt

CAS 103213-51-0

Couple Type Modified nucleotides

Formula $C_{12}H_{11}N_5O_6P \cdot Na$

MW 375.2

Purity > 98% HPLC

Quantity 10µmol≈3.8mg

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-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-de@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-uk@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com