

8-N₃-ε-cAMP

Immobilizable cAMP analogue for
fluorescent tagging

BLG-A077-05 (5 x 1 μmol pack size) is not sold in the U.S. or Canada.
Please [contact us](#) for available options.

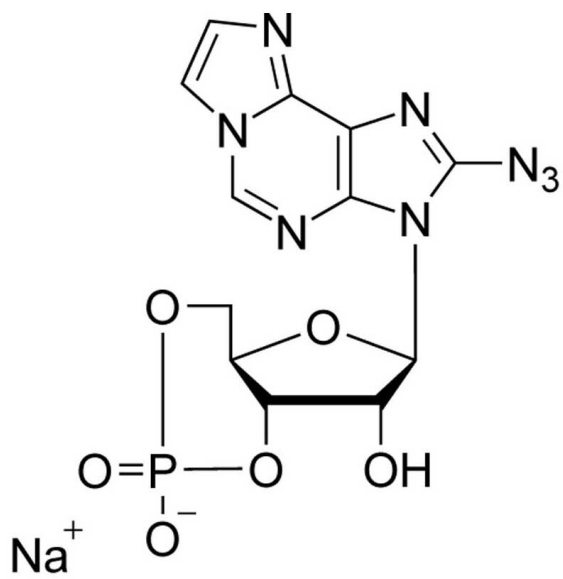
Ordering Information

[Order Online »](#)

BLG-A077-01	1 μmol
BLG-A077-05	5x1 μmol

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 8-Azido-1,N6-ethenoadenosine-3',5'-cyclic monophosphate . sodium salt

CAS 60902-12-7

Couple Type Modified nucleotides

Formula $C_{12}H_{10}N_8O_6P \cdot Na$

MW 416.2

Purity > 98% HPLC

Quantity 1µmol≈0.4mg

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](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)