

2'-ADOC-cAMP

cAMP analogue with a hydrophilic spacer for improved solubility in water, suitable as a ligand in affinity chromatography and for modification with fluorophores or other markers.

BLG-A146-25 (5 x 5 μ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

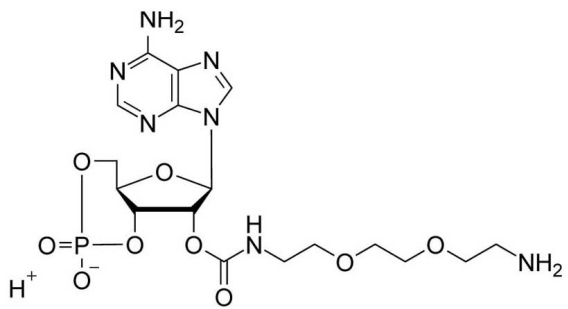
Ordering Information

[Order Online »](#)

| | |
|--------------------|---------------|
| BLG-A146-05 | 5 μ mol |
| BLG-A146-25 | 5x5 μ mol |

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 2'-O-(8-Amino-3,6-dioxaoctylcarbamoyl)adenosine-3',5'-cyclic monophosphate

Couple Type Modified nucleotides

Formula $C_{17}H_{26}N_7O_9P$

MW 503.4

Purity > 98% HPLC



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