

Sp-2-AEA-cAMPS

PDE-resistant Sp-cAMPS functionalized for immobilization or conjugation with markers and dyes

BLG-A093-25 (5 x 5 μ mol pack size) is not sold in the U.S. or Canada.

Please [contact us](#) for available options.

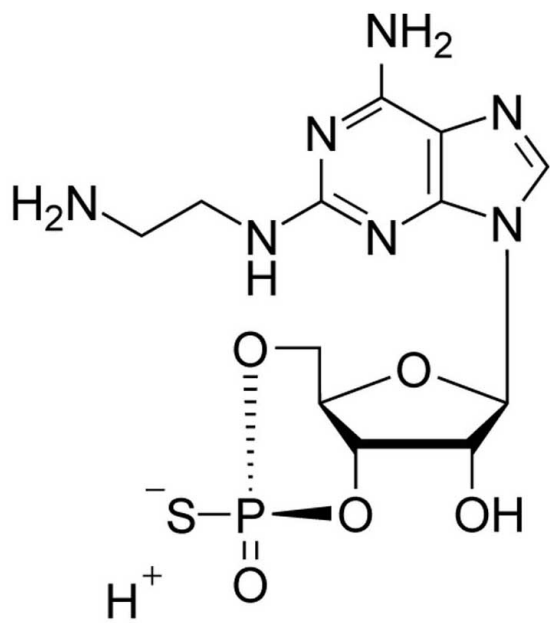
Ordering Information

[Order Online »](#)

| | |
|-------------|---------------|
| BLG-A093-05 | 5 μ mol |
| BLG-A093-25 | 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-(2-Aminoethylamino)adenosine-3',5'-cyclic monophosphorothioate, Sp-isomer

Couple Type Modified nucleotides

Formula $C_{12}H_{18}N_7O_5PS$

MW 403.4

Purity > 98% HPLC

Quantity 5µmol≈2mg



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