

Sp-2-Cl-cAMPS

Analogue of the protein kinase A activator Sp-cAMPS (Cat. No. A003) with increased lipophilicity and phosphodiesterase stability. The halogen in position 2 of the adenine nucleobase can be used for various substitutions including spacers and labels.

*BLG-C074-25 (5 x 5 μ mol pack size) is not sold in the U.S. or Canada.
Please [contact us](#) for available options.*

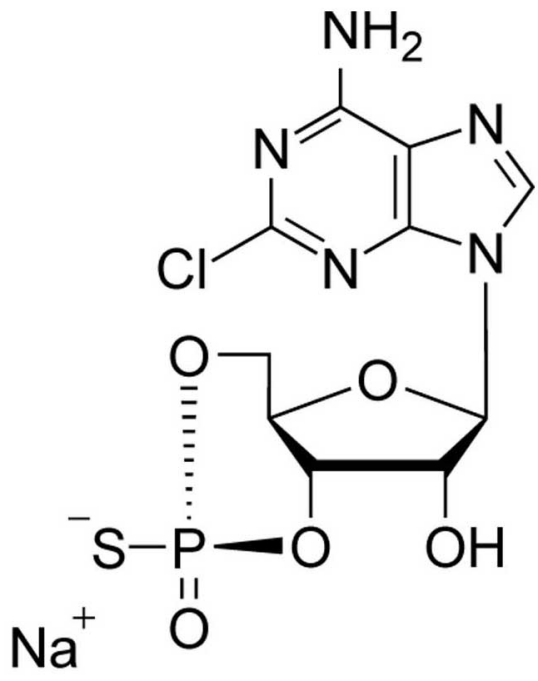
Ordering Information

[Order Online »](#)

BLG-C074-05	5 μ mol
BLG-C074-25	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-Chloroadenosine-3',5'-cyclic monophosphorothioate, Sp-isomer sodium salt

CAS 124854-63-3

Couple Type Modified nucleotides

Formula $C_{10}H_{10}ClN_5O_5PS \cdot Na$

MW 401.7

Purity > 99% 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