

Rp-c-diAMPS

Monophosphorothioate analogue of c-diAMP

Rp-c-diAMPS may be useful in studies on ligand-receptor interactions with c-diAMP and c-diGMP-binding proteins. The corresponding Sp-isomer Sp-c-diAMPS is also offered (Prod. No. BLG-C122).

BLG-C121-005 (5 x 0.1 μ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

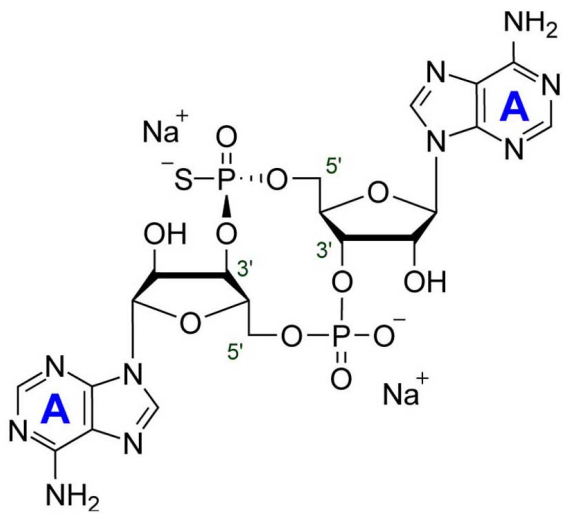
Ordering Information

[Order Online »](#)

BLG-C121-001	0.1 μ mol
BLG-C121-005	5x0.1 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 Cyclic diadenosine monophosphorothioate, Rp-isomer, sodium salt

Couple Type Modified nucleotide

Formula $C_{20}H_{24}N_{10}O_{11}P_2S$ (free acid)

MW 674.5 (free acid)

Purity > 95% HPLC

Quantity 0.1 μmol ≈ 67 μg

Solubility Soluble in water (≥3mM).

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-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