

Rp-cAMPS . triethyl ammonium salt

Membrane permeant and PDE-resistant inhibitor of PKA

Competitive inhibitor of protein kinase A type I and II (cyclic AMP antagonist). Resistant against mammalian cyclic nucleotide-dependent phosphodiesterases. Significantly more lipophilic compared to cyclic AMP.

BLG-A002T-25 (5 x 5 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

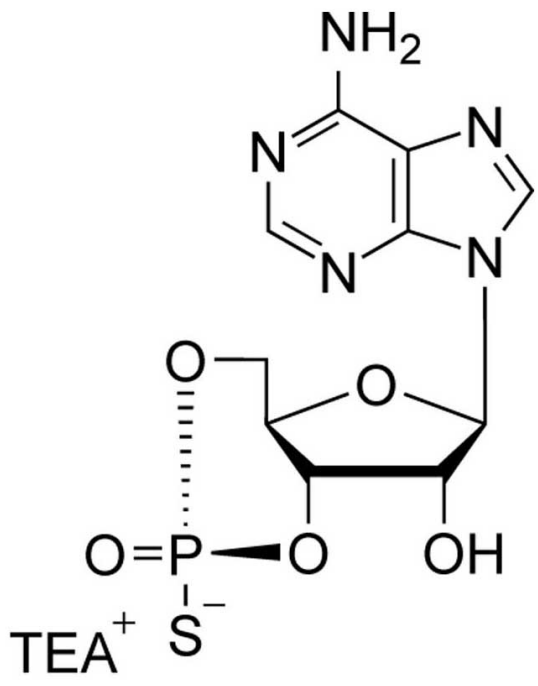
Ordering Information

[Order Online »](#)

BLG-A002T-05	5µmol
BLG-A002T-25	5x5µmol

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 Adenosine-3',5'-cyclic monophosphorothioate . triethyl ammonium salt, Rp-isomer

Appearance Lyophilized or crystallized sodium salt.

CAS 151837-09-1

Couple Type Inhibitor, Modified nucleotides

Formula $C_{10}H_{11}N_5O_5PS \cdot C_6H_{16}N$

MW 446.5

Purity > 98% HPLC

Quantity 5µmol≈2.2mg

Reconstitution Rinse tube walls carefully and preferably use ultrasonic or vortex to achieve total and uniform mixing.

Solubility Soluble in water or buffer.

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)