

Sp-cAMPS . sodium salt

PDE-resistant PKA agonist

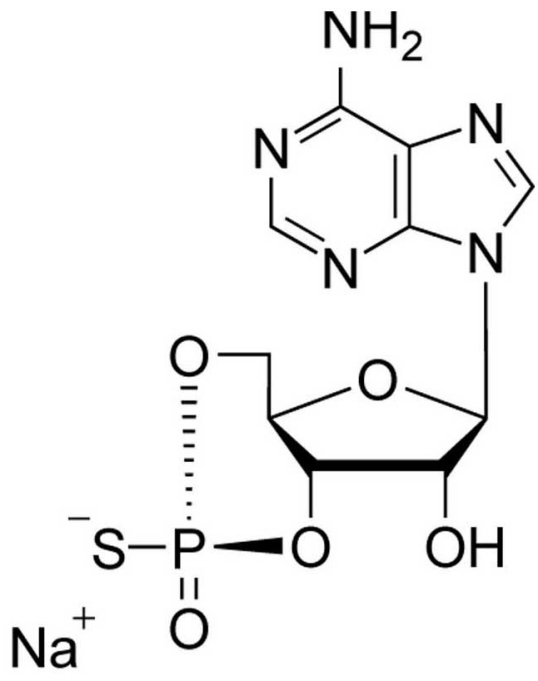
Activator of protein kinase A (cyclic AMP agonist). Considerable higher resistance against cyclic nucleotide phosphodiesterases compared to dibutyryl or 8-bromo cyclic AMP, no metabolic side effects. Membrane permeability comparable to 8-Br-cAMP.

BLG-A003S-50 (5 x 10 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

Ordering Information [Order Online »](#)

BLG-A003S-10	10µmol
BLG-A003S-50	5x10µmol

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 . sodium salt, Sp-isomer

CAS 71774-13-5

Couple Type Modified nucleotides

Formula $C_{10}H_{11}N_5O_5PS \cdot Na$

MW 367.3

Purity > 99% HPLC

Quantity 10µmol≈3.7mg

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