

The second messenger cyclic AMP activates protein kinase A, cAMP-gated ion channels, the CAP protein and some other target receptors. It is metabolized rapidly by cyclic nucleotide-dependent phosphodiesterases.

BLG-A001S-2500 (5 x 500 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

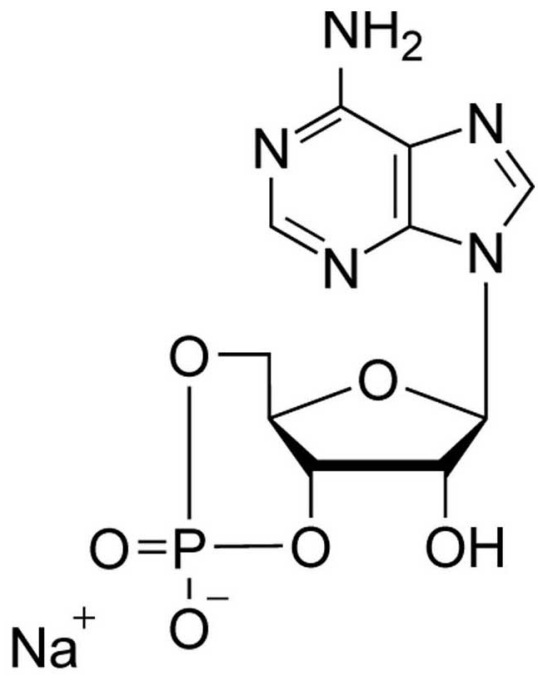
Ordering Information

[Order Online »](#)

| | |
|-----------------------|-----------|
| BLG-A001S-2500 | 5x500µmol |
|-----------------------|-----------|

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 monophosphate . sodium salt

CAS 37839-81-9

Couple Type Modified nucleotides

Formula $C_{10}H_{11}N_5O_6P \cdot Na$

MW 351.2

Purity > 99% HPLC

Quantity 500µmol≈176mg

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