

cAMP-AM

Metabolically activatable prodrug of the second messenger cyclic AMP

A membrane permeant precursor of the second messenger cyclic AMP. Releases cAMP after metabolic activation by esterases which is trapped inside the cell and metabolized quickly yielding a pulse-type signal.

BLG-A022-05 (5 x 1 μ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

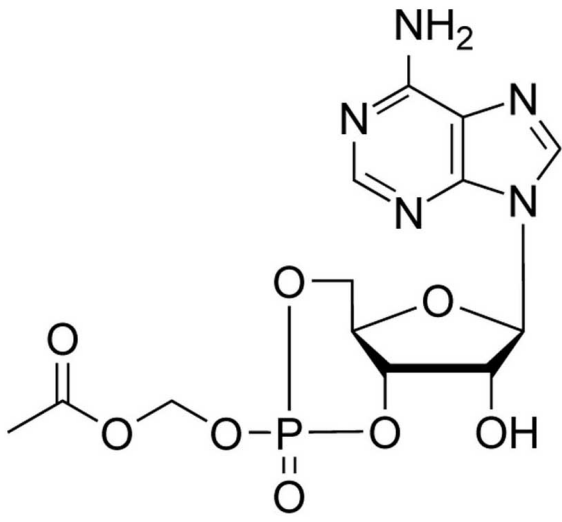
Ordering Information

[Order Online »](#)

BLG-A022-01	1 μ mol
BLG-A022-05	5x1 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 Adenosine-3',5'-cyclic monophosphate . acetoxymethyl ester

CAS 159764-93-9

Couple Type Modified nucleotides

Formula $C_{13}H_{16}N_5O_8P$

MW 401.3

Purity > 97% HPLC for mixture of isomers

Quantity 1µmol≈0.4mg

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