

8-ABA-cAMP

cyclic AMP functionalized for conjugation
with markers and dyes or for immobilization

Site selective cAMP analogue with preference for site B of R I of cAMP-dependent protein kinase. In combination with 8-piperidino-cAMP specific synergistic activation of cAK type I can be obtained. Also suitable as a ligand in affinity chromatography and for modification with fluorophores or other markers.

*BLG-A017-25 (5 x 5 μ mol pack size) is not sold in the U.S. or Canada.
Please [contact us](#) for available options.*

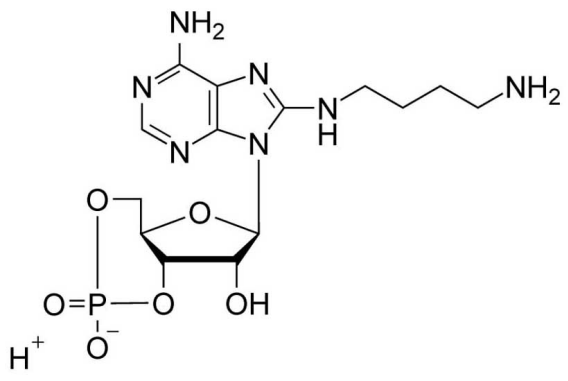
Ordering Information

[Order Online »](#)

BLG-A017-25	5x5 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 8-(4-Aminobutyl)aminoadenosine-3',5'-cyclic monophosphate

CAS 82927-68-2

Couple Type Modified nucleotides

Formula $C_{14}H_{22}N_7O_6P$

MW 415.4

Purity > 98% HPLC

Quantity 5µmol≈2.1mg

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