

2-ADOA-cAMP

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

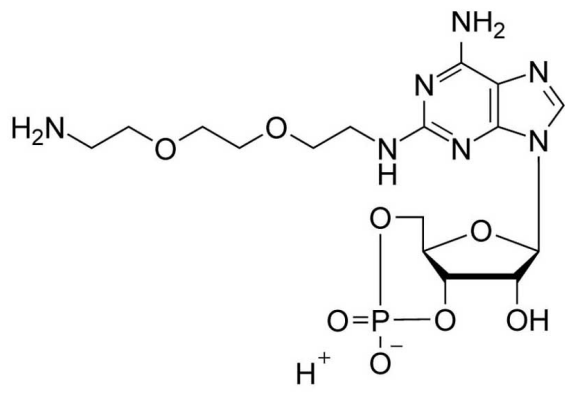
Due to its free aliphatic amino group various activated structures such as fluorescent dyes or other labels can be attached to 2-ADOA-cAMP. Another application is its use as a ligand for affinity chromatography of cAMP- and cGMP-dependent protein kinases and phosphodiesterases.

BLG-A136-25 (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-A136-05	5 μ mol
BLG-A136-25	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 2-(8-Amino-3,6-dioxaoctylamino)adenosine-3',5'-cyclic monophosphate

Couple Type Modified nucleotides

Formula $C_{16}H_{26}N_7O_8P$

MW 475.4

Purity > 98% HPLC

Quantity 5µmol≈2.4mg

Solubility Due to its ability to form internal and polymeric salts 2-ADOA-cAMP has limited solubility in water or buffer, however, a ≥13mM solution in water can be achieved. The compound is better soluble in dilute alkali of pH 9.5 and can, after dissolution, be titrated back to neutral. In addition, gentle heating and an ultrasonic bath usually helps to get complete dissolution.

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