

2-Biotin-15-cAMP

Cyclic AMP conjugated with a biotin label

2-Biotin-AHA-cAMP is an analogue of the natural signal molecule cyclic AMP in which the hydrogen in position 2 of the heterocyclic nucleobase is replaced by a long spacer, carrying a biotin at its end.

- Analogue of cyclic AMP with biotin conjugate most probably suitable as tracer in immunoassays for the measurement of cAMP concentrations,
- biotin as primary probe allows for flexible end point determinations with commercially available labeled streptavidin derivatives.

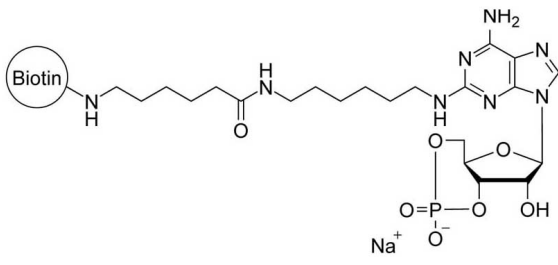
In spite of its modification, 2-Biotin-AHA-cAMP could be still sensitive against phosphodiesterases. For corresponding PDE-resistant structures, such as phosphorothioate-modified congeners or other types of spacers please inquire.

BLG-B041-05 (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-B041-01	1 μ mol
BLG-B041-05	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 2-(6-(6-[Biotinyl]aminohexanoyl)aminohexylamino)adenosine-3',5'-cyclic monophosphate . sodium salt

Appearance Crystallized or lyophilized salt.

Couple Type Modified nucleotides

Formula $C_{32}H_{50}N_{10}O_9PS \cdot Na$

MW 804.8

Purity > 98% HPLC

Quantity 1µmol≈0.8mg

Solubility Soluble in water.

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