

5'-Biotin-16-2',3'-cAMP

Biotin conjugated to 2',3'-cAMP

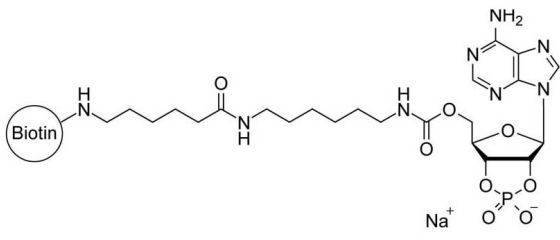
Biotinylated analogue of the signalling nucleotide 2',3'-cAMP (Biolog Cat. No. A 307) in which the biotin label is attached to the ribose 5' position via a 9-atom spacer. Potential tracer in immunoassays, for affinity chromatography or pull-down assays.

*BLG-B267-05 (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-B267-01	1 μ mol
BLG-B267-05	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 5'-O-(6-(6-[Biotinyl]aminohexanoyl)aminohexylcarbamoyl)adenosine-2',3'-cyclic monophosphate . sodium salt, 5'-[Biotin]-AHC-2',3'-cAMP

Couple Type Modified nucleotides

Formula $C_{33}H_{50}N_{10}O_{10}PS$

MW 832.85

Purity > 97% HPLC

Quantity 1µmol is equivalent to ~0.8 mg

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