

# 4-Biotin-14-cCMP

Cyclic CMP conjugated with a biotin label

The biotin label is connected to the N<sup>4</sup>-position of cyclic CMP (cCMP, Cat. No. C 001) via a 14-atom spacer. 4-Biotin-14-cCMP may be useful as tracer in immunoassays, for affinity chromatography or pull-down assays.

*BLG-B220-05 (5 x 1 µmol pack size) is not sold in the U.S. or Canada.  
Please [contact us](#) for available options.*

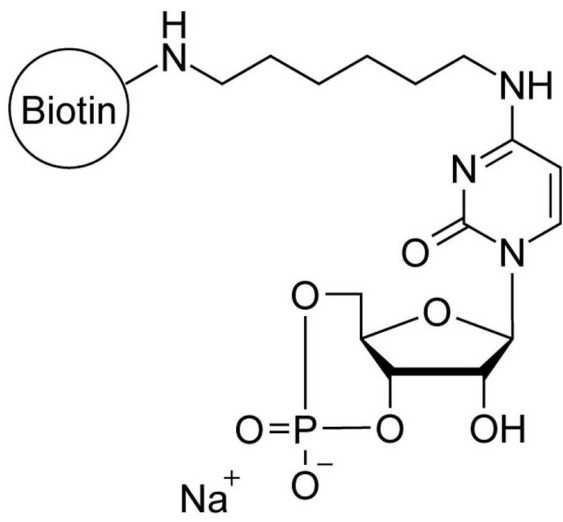
## Ordering Information

[Order Online »](#)

BLG-B220-01	1µmol
BLG-B220-05	5x1µmol

## 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** N4-(6-(6-[Biotinyl]aminohexanoyl)aminohexyl)cytidine-3',5'-cyclic monophosphate . sodium salt, 4-[Biotin]-AH-cCMP

**Appearance** Solid.

**Couple Type** Modified nucleotides

**Formula**  $C_{31}H_{49}N_7O_{10}PS \cdot Na$

**MW** 765.8

**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](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)