

8-N -6-Biotin-10-cAMP

Analogue for photoaffinity labelling of cAMP binding proteins

BLG-A127-005 (5 x 0.1 μ mol pack size) is not sold in the U.S. or Canada.
Please [contact us](#) for available options.

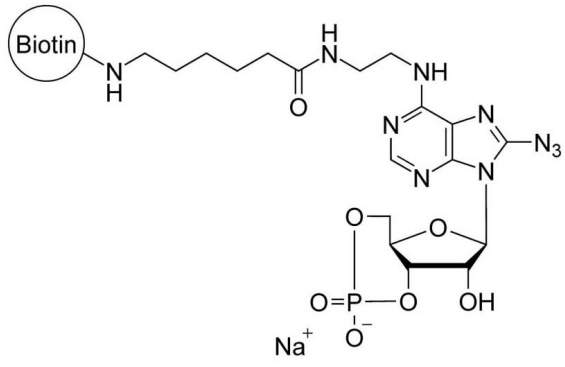
Ordering Information

[Order Online »](#)

BLG-A127-001	0.1 μ mol
BLG-A127-005	5x0.1 μ mol

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	8-Azido-N6-(2-(6-[biotinyl]amino)hexanoyl)aminoethyl)adenosine-3',5'-cyclic monophosphate . sodium salt
Couple Type	Modified nucleotides
Formula	$C_{28}H_{12}NO_9PS \cdot Na$
MW	774.7
Purity	> 98%
Quantity	0.1 μ mol \approx 80 μ g
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