

2',2''-Di-Biotin-16-c-diAMP

Dibiotinylated analogue of the bacterial second messenger c-diAMP

2',2''-Di-[Biotin]-AHC-c-diAMP is an analogue of the bacterial second messenger c-diAMP (BIOLOG Cat. No. BLG-C088) in which biotin moieties have been attached to both ribose 2'-hydroxy groups via 16-atom spacers.

BLG-D122-005 (5 x 0.1 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

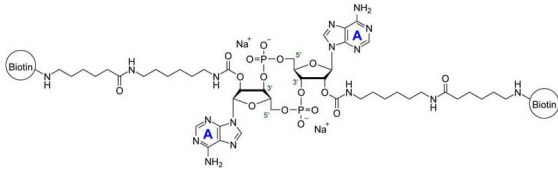
Ordering Information

[Order Online »](#)

BLG-D122-001	0.1µmol
BLG-D122-005	5x0.1µmol

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', 2''-O-(Di-(6-(6-[biotinyl]aminohexanoyl)aminohexylcarbamoyl))-cyclic diadenosine monophosphate . sodium salt

Couple Type Modified nucleotides

Formula $C_{66}H_{102}N_{20}O_{20}P_2S_2$ (free acid)

MW 1621.7 (free acid)

Purity > 95% HPLC

Quantity 0.1 µmol ≈ 0.2mg

Solubility Soluble in water or in aqueous buffers (>1mM).

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