

2',2''-Di-AHC-c-diGMP

Functionalized c-diGMP for labelling or immobilization

Analogue of the bacterial second messenger c-diGMP (Cat. No. BLG-C057) which is suitable as a ligand in affinity chromatography or for coupling of various labelling structures including fluorophores.

BLG-D112-005 (5 x 0.1 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

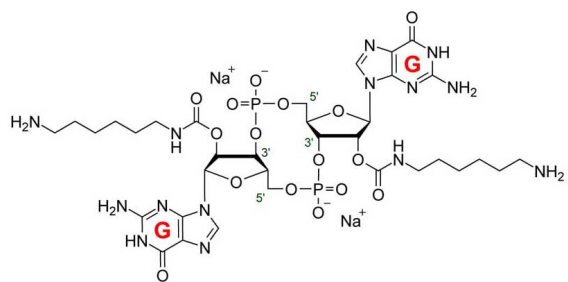
Ordering Information

[Order Online »](#)

BLG-D112-001	0.1µmol
BLG-D112-005	5x0.1µmol

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	2',2"-Di-AHC-c-diAMP has sufficient stability at room temperature and does not need special care during handling or shipment. Nevertheless, we recommend that the compound should be stored in the freezer, for longer storage periods preferably in freeze-dried form.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	2', 2"-O-(Di-[6-aminohexylcarbamoyl])-cyclic diguanosine monophosphate . sodium salt
Couple Type	Modified nucleotides
Formula	$C_{34}H_{52}N_{14}O_{16}P_2$ (free acid)
MW	974.8 (free acid)
Purity	> 95% HPLC
Quantity	0.1 μmol ≈ ~97 μg
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-fr@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