

2'-AHC-c-diGMP

Functionalized c-diGMP for labelling or immobilization

Precursor for modification with fluorophores and other markers. Also suitable as a ligand for immobilization to yield affinity gels.

BLG-A151-005 (5 x 0.1 μ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

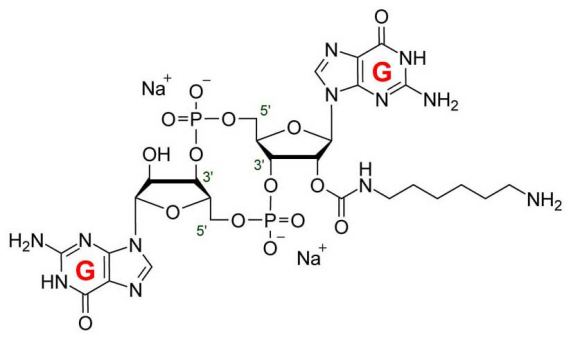
Ordering Information

[Order Online »](#)

BLG-A151-001	0.1 μ mol
BLG-A151-005	5x0.1 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 2'-O-(6-Aminohexylcarbamoyl)cyclic diguanosine monophosphate . sodium salt

CAS 1422761-54-3

Couple Type Modified nucleotides

Formula $C_{27}H_{38}N_{12}O_{15}P_2$

MW 832.6 (free acid)

Purity > 95% HPLC

Quantity 0.1 μmol ≈ 83 μg

Solubility Soluble in water or aqueous buffers.

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