

2'-[sCya3]-AHC-c-diGMP

Fluorescent analogue of c-diGMP

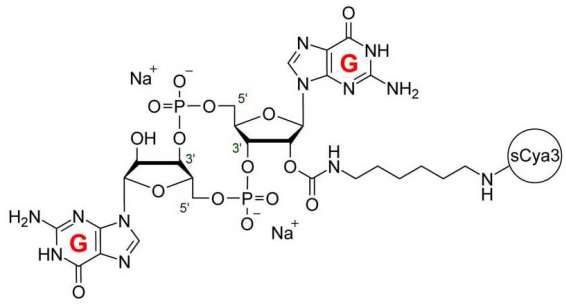
2'-[sCya3]-AHC-c-diGMP is a fluorescently labelled version of the bacterial cyclic dinucleotide second messenger c-diGMP (Cat. No. C 057), fluorescent with λ_{exc} 548 nm and λ_{em} 563 nm. The sulfo-Cyanine 3 dye is connected to one of the ribose-2' positions of c-diGMP via a 9 atom spacer.

BLG-C412-005 (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-C412-001	0.1 μ mol
BLG-C412-005	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-[sulfo-Cyanine3]-aminohexylcarbamoyl)-cyclic diguanosine monophosphate . sodium salt

Appearance Solid

Couple Type Modified nucleotides

Formula $C_{57}H_{72}N_{14}O_{22}P_2S_2$ (free acid)

MW 1431.4

Purity > 95% HPLC

Quantity 0.1 µmol is equivalent to ~0.1 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

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