

c[3'-[sCya5]-AHC-G(2',5')pA(3',5')p]

Fluorescently labelled 2'3'-cGAMP

Fluorescently labelled version of the metazoan cyclic dinucleotide second messenger and STING activator c[G(2',5')pA(3',5')p] (aka cGAMP(2'-5') or 2'3'-cGAMP, Cat. No. C 161), fluorescent with λ_{exc} 644 nm and λ_{em} 664 nm. c[G(2',5')pA(3',5')p] was found to be the metazoan second messenger produced by the mammalian innate immune DNA sensor cGAMP synthase (cGAS).

BLG-C270-005 (5 x 0.1 μ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

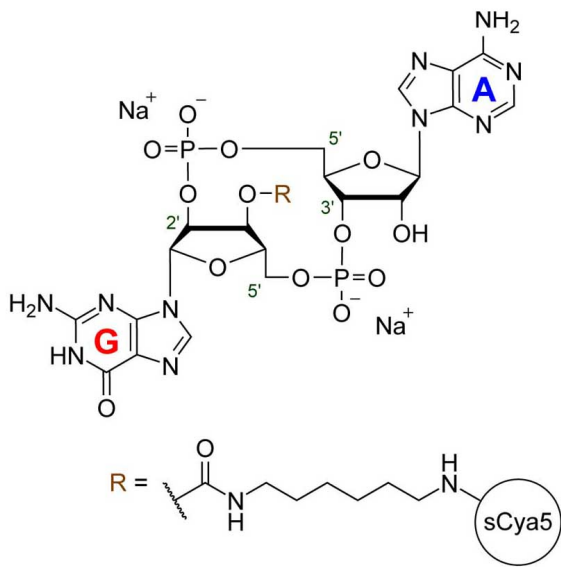
Ordering Information

[Order Online »](#)

BLG-C270-001	0.1 μ mol
BLG-C270-005	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 Cyclic (3'-O-(6-[sulfo-cyanine5]amino-hexyl-carbamoyl)guanosine-(2'→5')-monophosphate-adenosine-(3'→5')-monophosphate) . sodium salt

Couple Type Modified nucleotides

Formula $C_{59}H_{74}N_{14}O_{21}P_2S_2$ (free acid)

MW 1441.4 (free acid)

Purity > 95% HPLC

Quantity 0.1 μmol is equivalent to ~144 μg



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