

c[3'-[sCya7]-AHC-G(2',5')pA(3',5')p]

Fluorescent analogue of 2',3'-cGAMP

BLG-C240-005 (5 x 0.1 μ mol pack size) is not sold in the U.S. or Canada.

Please [contact us](#) for available options.

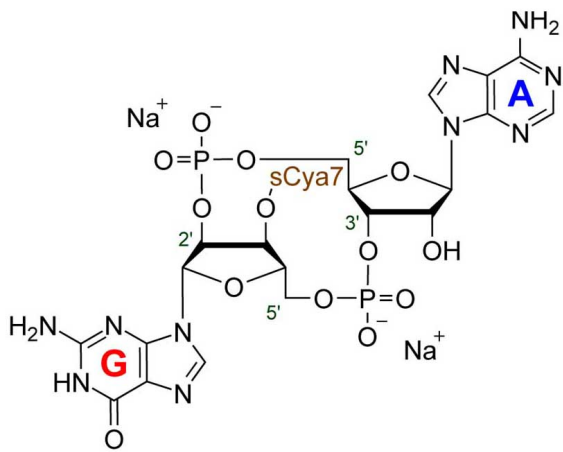
Ordering Information

[Order Online »](#)

BLG-C240-001	0.1 μ mol
BLG-C240-005	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-cyanine7]amino)hexylcarbamoyl)guanosine-(2'→5')-monophosphate-adenosine-(3'→5')-monophosphate, sodium salt

Formula $C_{64}H_{80}N_{14}O_{21}P_2S_2$

MW 1507.5

Purity > 95% HPLC



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