

c[G(2',5')pS-A(3',5')pS], isomer 1 and isomer 2

Set of two isomeric di-thiophosphate analogues of the metazoan cyclic

dinucleotide second messenger 2'3'-cGAMP
This set contains two isomeric di-thiophosphate analogues of the metazoan cyclic dinucleotide second messenger c[G(2',5')pA(3',5')p] (2'3'-cGAMP, Cat. No. BLG-C161).

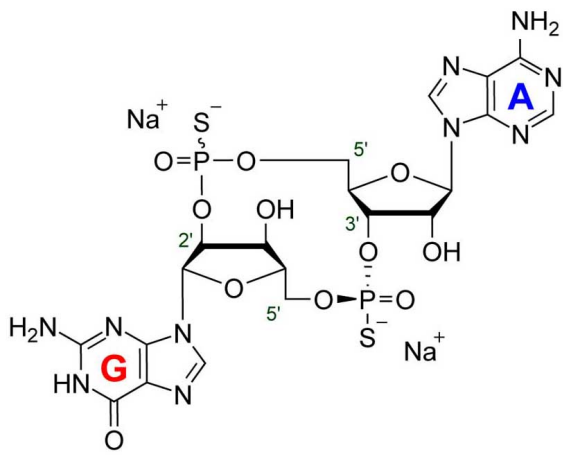
Ordering Information

[Order Online »](#)

BLG-SET101	2x0.1µmol
------------	-----------

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(guanosine-(2'→5')-monophosphorothioate-adenosine-(3'→5')-monophosphorothioate), 2'3'-cGAMPSS, 2'3'-cGsAsMP

CAS 1638242-56-4/1638243-00-1

Couple Type Modified nucleotides

Formula $C_{20}H_{24}N_{10}O_{11}P_2S_2$

Purity > 95% HPLC

Quantity 2×0.1µmol≈71µg

Solubility Each isomer of c[G(2',5')pS-A(3',5')pS] is soluble to ≥ 1.4 mM in water. Please rinse tube walls carefully and preferably use ultrasonic or vortex to achieve total and uniform mixing. When opening the tube please make sure that no substance is lost within the cap.

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