

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

BLG-C247-005 (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 »](#)

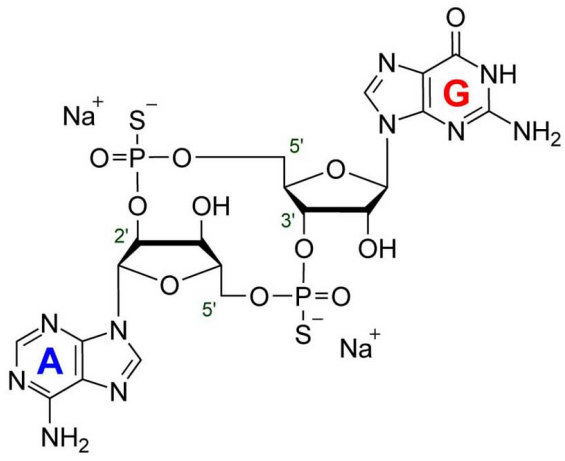
BLG-C247-005

5x0.1 μ mol

Manuals, SDS & CofA

[View Online »](#)

c[A(2',5')pS-G(3',5')pS], isomer 2 (syn. 3'2'-cGAMPSS, isomer 2) is an isomeric di-thiophosphate derivative of the recently discovered nucleotide second messenger c[A(2',5')pG(3',5')p] (3'2'-cGAMP, Cat. No. C 238). The di-thiophosphate modifications are expected to render the compound resistant to degradation by 3'2'-cGAMP hydrolysing enzymes. c[A(2',5')pS-G(3',5')pS], isomer 1 is available as well (syn. 3'2'-cGAMPSS, isomer 1, Cat. No. C 246).



Handling & Storage

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 3'2'-cGAMPSS, isomer 2, 3',5'-2',5'-cGAMPSS, Cyclic (adenosine-(2'→5')-monophosphorothioate-guanosine- (3'→5'-monophosphorothioate), isomer 2 . sodium salt

Excitation Maximum 256 nm

Formula $C_{20}H_{24}N_{10}O_{11}P_2S_2$ (free acid)

MW 706.6 (free acid)

Purity > 95% HPLC

Quantity 0.1 µmol / ~ 71 µg

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