

$[^{13}\text{C}_{20}, ^{15}\text{N}_{10}]$ - c[G(2',5')pA(3',5')p] / [$^{13}\text{C}_{20}, ^{15}\text{N}_{10}$]-2'3'- cGAMP

BLG-C410-0025 (5 x 0.05 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

Stable isotope version of the metazoan cyclic dinucleotide second messenger

c[G(2',5')pA(3',5')p] / 2'3'-cGAMP

Ordering Information

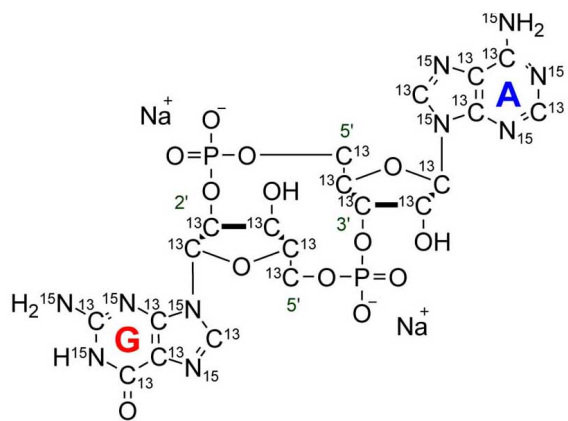
[Order Online »](#)

BLG-C410-0005	0.05µmol
BLG-C410-0025	5x0.05µmol

Manuals, SDS & CofA

[View Online »](#)

$[^{13}\text{C}_{20}, ^{15}\text{N}_{10}]$ -c[G(2',5')pA(3',5')p] is the stable isotope version of the metazoan cyclic dinucleotide second messenger c[G(2',5')pA(3',5')p] / 2'3'-cGAMP (Biolog Cat. No. C 161). c[G(2',5')pA(3',5')p] is produced by the mammalian innate immune DNA sensor cGAMP synthase (cGAS) and potentially activates the immune signalling protein STING (stimulator of interferon genes). $[^{13}\text{C}_{20}, ^{15}\text{N}_{10}]$ -c[G(2',5')pA(3',5')p] is suitable as internal standard in LC-MS applications for c[G(2',5')pA(3',5')p]. The isotopic purity is 98% for ^{13}C and ^{15}N , respectively.



Handling & Storage

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Cyclic (guanosine-[¹³C₁₀,¹⁵N₅]-[2'→5']-monophosphate-adenosine- [¹³C₁₀,¹⁵N₅]-[3'→5']-monophosphate) . sodium salt, [¹³C₂₀,¹⁵N₁₀]-cGAMP(2'-5'), [¹³C₂₀,¹⁵N₁₀]-2',5'-3',5'-cGAMP

Formula $^{13}\text{C}_{20}\text{H}_{24}^{15}\text{N}_{10}\text{O}_{13}\text{P}_2$ (free acid)

MW 704.4 (free acid)

Purity > 95% HPLC

Quantity 0.05µmol = ~35µg

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