

[¹³C₂₀, ¹⁵N₁₀]-c-diGMP

Stable isotope version of the bacterial second messenger c-diGMP

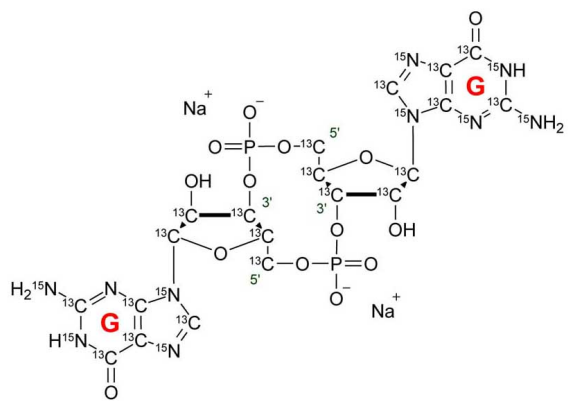
Stable isotope version of the bacterial second messenger and STING ligand c-diGMP (Cat. No. BLG-C057) with 98% isotopic purity of ¹³C and ¹⁵N, respectively. Suitable as internal standard in LC-MS applications for c-diGMP.

BLG-C183-0025 (5 x 0.05 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

Ordering Information [Order Online »](#)

BLG-C183-0005	0.05µmol
BLG-C183-0025	5x0.05µmol

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 diguanosine-[13C20,15N10]-monophosphate . sodium salt

Appearance Solid

Formula $^{13}\text{C}_{20}\text{H}_{24}^{15}\text{N}_{10}\text{O}_{14}\text{P}_2$ (free acid)

MW 720.4 (free acid)

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