

[¹⁵N₁₀]-c-diAMP

Stable isotope version of the bacterial second messenger c-diAMP

BLG-C409-0025 (5 x 0.05 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

Ordering Information

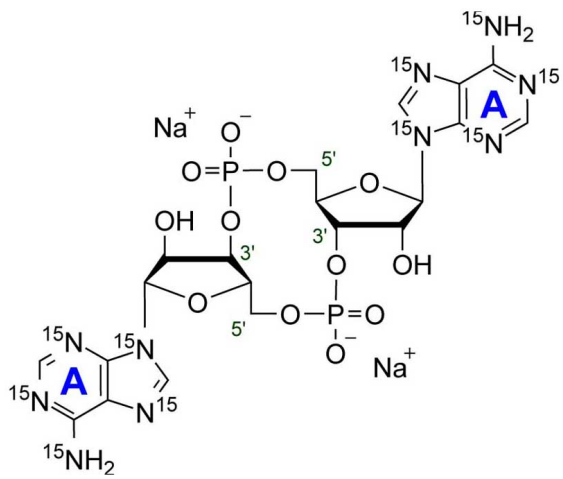
[Order Online »](#)

BLG-C409-0005	0.05µmol
BLG-C409-0025	5x0.05µmol

Manuals, SDS & CofA

[View Online »](#)

[¹⁵N₁₀]-c-diAMP is the stable isotope version of the bacterial second messenger c-diAMP (Biolog Cat. No. C 088) that was found to be involved in a wide variety of physiological processes in bacteria such as biofilm formation, maintenance of osmotic pressure and acid stress resistance. [¹⁵N₁₀]-c-diAMP is suitable as internal standard in LC-MS applications for c-diAMP. The isotopic purity is 98%.



Handling & Storage

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Cyclic diadenosine-[15N10]-monophosphate . sodium salt

Formula $C_{20}H_{24}^{15}N_{10}O_{12}P_2$ (free acid)

MW 668.4 (free acid)

Purity > 95% HPLC

Quantity 0.05µmol = ~33µg

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