

c-diAMP

Bacterial second messenger

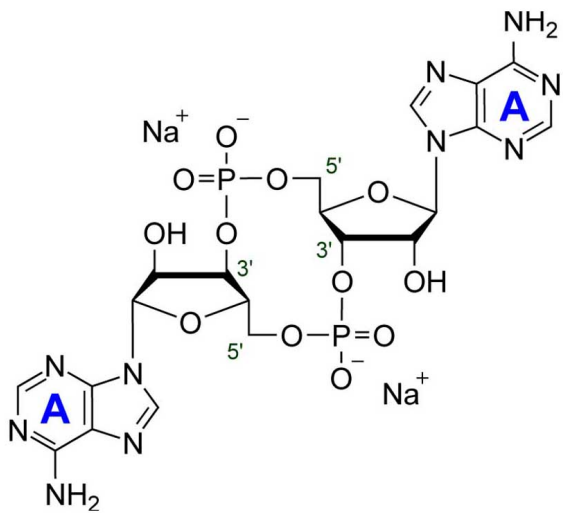
c-diAMP may be an additional purine-based signalling nucleotide in prokaryotes. The structurally related bacterial second messenger cyclic diguanosine monophosphate (c-diGMP) is offered as well (Prod. No. [BLG-C057](#)).

*BLG-C088-05 (5 x 1 µmol pack size) is not sold in the U.S. or Canada.
Please [contact us](#) for available options.*

Ordering Information [Order Online »](#)

BLG-C088-01	1 µmol
BLG-C088-05	5x1 µmol

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 diadenosine monophosphate, Cyclic bis (3'→5') diadenylic acid . disodium salt

CAS 54447-84-6

Couple Type Modified nucleotides

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

MW 658.4 (free acid)

Purity > 98% HPLC (typically > 99%)

Quantity 1 µmol = ~0.7 mg

Solubility Soluble in water (≥5mM).

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