

c-(ϵ -ApGp)

BLG-C226-005 (5 x 0.1 μ mol pack size) is not sold in the U.S. or Canada.
Please [contact us](#) for available options.

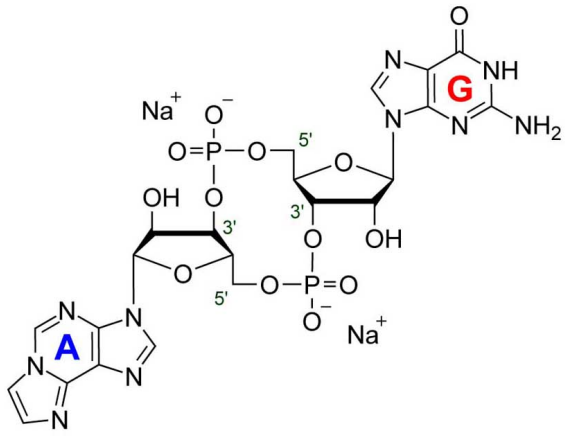
Ordering Information

[Order Online »](#)

BLG-C226-001	0.1 μ mol
BLG-C226-005	5x0.1 μ mol

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Cyclic (1, N6-ethenoadenosine monophosphate-guanosine monophosphate) . sodium salt, ϵ -3'-3'-cGAMP, c[ϵ -A(3',5')pG(3',5')p], c[G(3',5')p- ϵ -A(3',5')p]
Concentration	0.1 μ mol= \sim 70 μ g
Formula	$C_{22}H_{24}N_{10}O_{13}P_2$ (free acid)
MW	698.4
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