

c-tetraGMP

c-tetraGMP is a cyclic nucleotide in which four 5'-GMP units are interconnected via 3'-5' phosphodiester bonds to form a cyclic structure. c-tetraGMP is structurally related to the novel bacterial second messenger c-tetraAMP (Cat. No. C 335) which is involved in the Type III CRISPR-Cas-associated detection and degradation of invasive genetic elements in many prokaryotes.

*BLG-C359-025 (5 x 0.5 μ mol pack size) is not sold in the U.S. or Canada.
Please [contact us](#) for available options.*

Ordering Information

[Order Online »](#)

BLG-C359-005	0.5 μ mol
BLG-C359-025	5x0.5 μ mol

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 tetraguanosine monophosphate . sodium salt

Couple Type Modified nucleotides

Formula $C_{40}H_{48}N_{20}O_{28}P_4$ (free acid)

MW 1380.8 (free acid)

Purity > 95% HPLC

Quantity 0.5µmol is equivalent to ~0.7 mg



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