

c[3'-MANT- G(2',5')pA(3',5')p]

Fluorescent analogue of 2',3'-cGAMP

Fluorescent analogue of the metazoan cyclic dinucleotide second messenger and STING activator c[G(2',5')pA(3',5')p] (aka cGAMP(2'-5') or 2'3'-cGAMP, Cat. No. BLG-C161). The MANT fluorophore (λ_{exc} 355 nm; λ_{em} 448 nm) has a certain sensitivity for its environment and can change its spectral properties upon binding.

BLG-C300-005 (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-C300-001 | 0.1 μ mol |
| BLG-C300-005 | 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 (3'-O-(N'- methylanthraniloyl)guanosine-(2'→5')-monophosphate-adenosine-(3'→5')-monophosphate . sodium salt |
| Appearance | Solid |
| Formula | $C_{28}H_{31}N_{11}O_{14}P_2$ (free acid) |
| MW | 807.6 (free acid) |
| Purity | > 95% HPLC |
| 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