

c-(Ap-8-Fluo-AET-Gp)

Fluorescent analogue of cGAMP

*BLG-C159-005 (5 x 0.1 μ mol pack size) is not sold in the U.S. or Canada.
Please [contact us](#) for available options.*

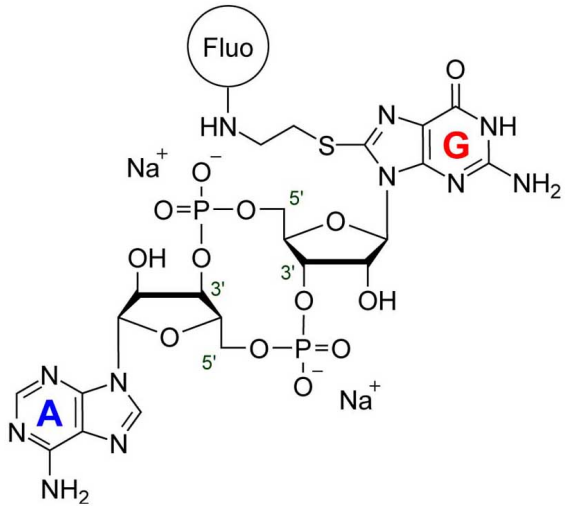
Ordering Information

[Order Online »](#)

BLG-C159-001	0.1 μ mol
BLG-C159-005	5x0.1 μ mol

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	Light-sensitive
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Cyclic (adenosine monophosphate-8-(2-[fluoresceinyl]aminoethylthio)guanosine monophosphate) . sodium salt
Formula	$C_{43}H_{39}N_{11}O_{19}P_2S$ (free acid)
MW	1107.9 (free acid)
Purity	> 95% HPLC
Quantity	0.1 μ mol \approx ~0.1 mg
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