

c[8-AET- G(2',5')pA(3',5')p]- Agarose

2',5'-cGAMP immobilized on agarose gel

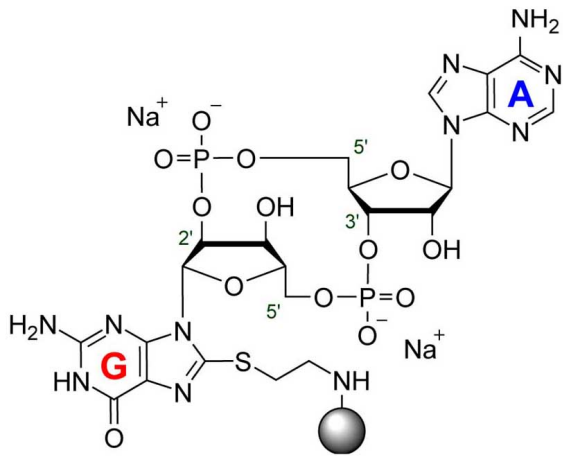
Ordering Information

[Order Online »](#)

BLG-C184-06	0.6ml
BLG-C184-25	2.5ml
BLG-C184-60	6ml

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	Stable at room temp but can be kept in freezer for long term storage in freeze dried form
Long Term Storage	+4°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Cyclic (8-(2-aminoethylthio)guanosine-(2'→5')-monophosphate-adenosine-(3'→5')-monophosphate), immobilized on agarose gel
Couple Type	Modified nucleotides
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