

Rp-8-AET-cGMPS-Agarose

Immobilized cGMP antagonist

The metabolically stable protein kinase G inhibitor Rp-cGMPS immobilized on agarose by an aminoethylthio spacer attached to position 8 of the ligand. Its reduced binding affinity towards protein kinase G compared to cGMP should result in milder desorption conditions in affinity chromatography. The gel can be used for affinity chromatography of various cyclic nucleotide-responsive proteins such as protein kinases, phosphodiesterases and others. Gels with the corresponding Sp-isomer (Sp-8-AET-cGMPS-Agarose (Prod. No. [BLG-A010](#))) and the parent cyclic phosphate (8-AET-cGMP-Agarose (Prod. No. [BLG-A019](#))) are also available.

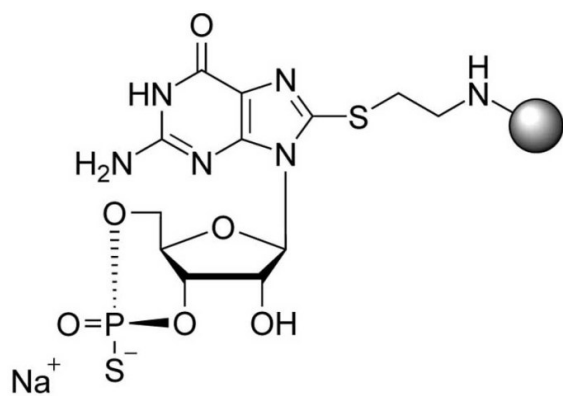
Ordering Information

[Order Online »](#)

BLG-A009-06	0.6ml
BLG-A009-25	2.5ml
BLG-A009-60	6ml

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Long Term Storage +4°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 8-(2-Aminoethyl)thioguanosine-3',5'-cyclic monophosphorothioate, Rp-isomer; immobilized on agarose gel

Couple Type Modified nucleotides

Formulation Liquid. 0.05M Na₂HPO₄ buffer (pH7), containing 0.1% sodium azide.

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](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)