

Rp-2'-AHC-cGMPS- Agarose

PDE-resistant Rp-cGMPS immobilized on agarose gel

The PDE-stable protein kinase G (PKG) inhibitor Rp-cGMPS immobilized on agarose by an aminohexylcarbamoyl spacer attached to the ribose 2' position of the ligand. The gel can be used for affinity chromatography of various cyclic nucleotide-responsive proteins, especially those which tolerate modification of the 2' ribose hydroxyl group, such as some phosphodiesterases. This type of gel is also available with immobilized PDE-resistant Sp-cGMPS (Sp-2'-AHC-cGMPS-Agarose (Prod. No. [BLG-A052](#))) and with normal cGMP (2'-AHC-cGMP-Agarose (Prod. No. [BLG-A059](#))).

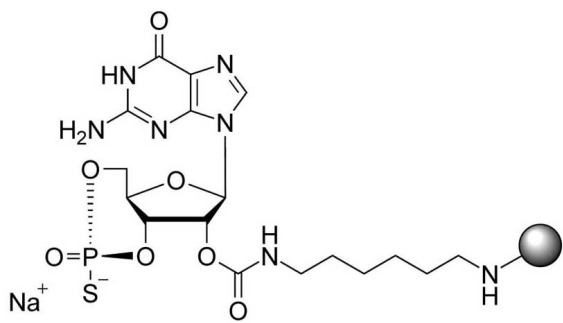
Ordering Information

[Order Online »](#)

BLG-A051-06	0.6ml
BLG-A051-25	2.5ml
BLG-A051-60	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 2'-(6-Aminohexylcarbamoyl)guanosine-3',5'-cyclic monophosphorothioate, Rp-isomer, 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