

Rp-8-AHA-cAMPS- Agarose

cyclic AMP antagonist immobilized on agarose gel

The metabolically stable protein kinase A inhibitor Rp-cAMPS immobilized on agarose by an aminohexylamino spacer attached to position 8 of the ligand. Free Rp-cAMPS binds to the protein kinase A holoenzyme but prevents dissociation to free catalytic and regulatory subunits. In addition, it is not hydrolyzed by mammalian phosphodiesterases. The gel can be used for affinity chromatography of various cyclic nucleotide-responsive proteins such as protein kinase A (presumably as its holoenzyme), phosphodiesterases and others. This ligand is also available with a shorter spacer arm (Rp-8-AEA-cAMPS-Agarose (Prod. No. [BLG-A007](#))).

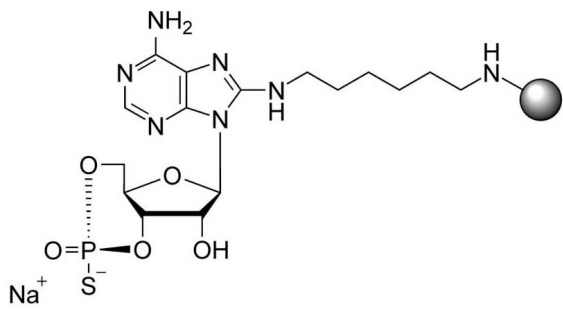
Ordering Information

[Order Online »](#)

BLG-A012-06	0.6ml
BLG-A012-25	2.5ml
BLG-A012-60	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-(6-Aminohexyl)aminoadenosine-3',5'-cyclic monophosphorothioate, Rp-isomer; immobilized on agarose gel

Couple Type Modified nucleotides

Formulation Liquid. 0.05M Na₂HPO₄ 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

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