

Rp-2'-AEC-cAMPS- Agarose

The protein kinase A inhibitor Rp-cAMPS (Prod. No. BLG-A002) immobilized on agarose by an aminoethylcarbamoyl spacer attached to the ribose 2'-position of the ligand. Free Rp-cAMPS is not metabolized by mammalian phosphodiesterases and thus represents an extraordinary stable ligand. In addition, its reduced binding affinity towards protein kinase A compared to cAMP results in milder desorption conditions during affinity chromatography. 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 offered with a longer spacer (Rp-2'-AHC-cAMPS-Agarose, Prod.. No. BLG-A049) as well

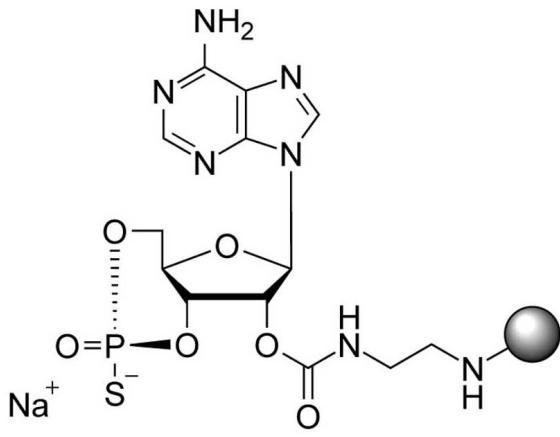
Ordering Information

[Order Online »](#)

BLG-A126-06	0.6ml
BLG-A126-25	2.5ml
BLG-A126-60	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'-O-(2-Aminoethylcarbamoyl)adenosine-3',5'-cyclic monophosphorothioate, Rp-isomer, immobilized on agarose gel

Couple Type Modified nucleotides



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