

Rp-8-AHA-cAMPS

The PKA inhibitor Rp-cAMPS functionalized for immobilization or conjugation with markers and dyes

The free terminal primary amino group in Rp-8-AHA-cAMPS, separated from the nucleotide by a hexyl spacer, is suitable for coupling to gels for affinity chromatography and for binding of various labels, e.g. fluorescent dyes. Rp-8-AHA-cAMPS is also available as ligand immobilized to agarose (Rp-8-AHA-cAMPS-Agarose, Prod. No. BLG-A012).

BLG-A085-25 (5 x 5 μ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

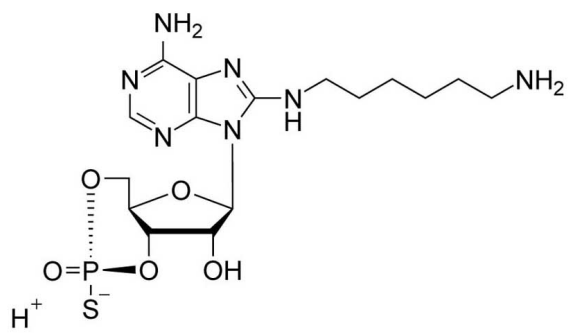
Ordering Information

[Order Online »](#)

BLG-A085-05	5 μ mol
BLG-A085-25	5x5 μ mol

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 8-(6-Aminohexylamino)adenosine-3',5'-cyclic monophosphorothioate, Rp-isomer

CAS 214272-02-3

Couple Type Modified nucleotides

Formula $C_{16}H_{26}N_7O_5PS$

MW 459.5

Purity > 98% HPLC

Quantity 5µmol≈2.3mg

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