

8-AHA-2'-O-Me-cAMP- Agarose

8-AHA-2'-O-Me-cAMP immobilized on agarose gel

Suitable for affinity chromatography of various cAMP-responsive proteins, especially those which tolerate modification of the ribose 2'-hydroxy group, such as the exchange protein activated by cyclic AMP (Epac) and certain phosphodiesterases.

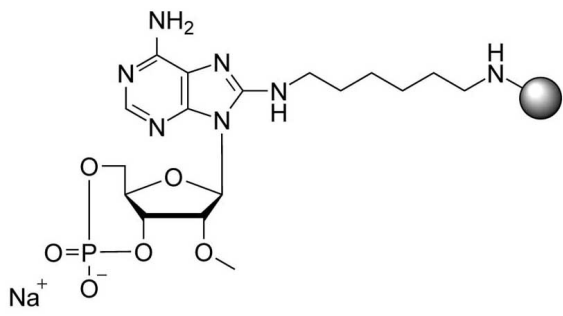
Ordering Information

[Order Online »](#)

BLG-A057-06	0.6ml
BLG-A057-25	2.5ml
BLG-A057-60	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)amino-2'-O-methyl-cAMP-Agarose

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