

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

The second messenger cyclic AMP with a methylated ribose 2'-hydroxyl group immobilized on agarose by an aminoethylthio spacer attached to position 8 of the ligand. Considered to be suitable for affinity chromatography of various cAMP-responsive proteins, especially those which tolerate modification of the ribose 2'-hydroxyl group, such as the exchange protein activated by cyclic AMP (Epac) and certain phosphodiesterases.

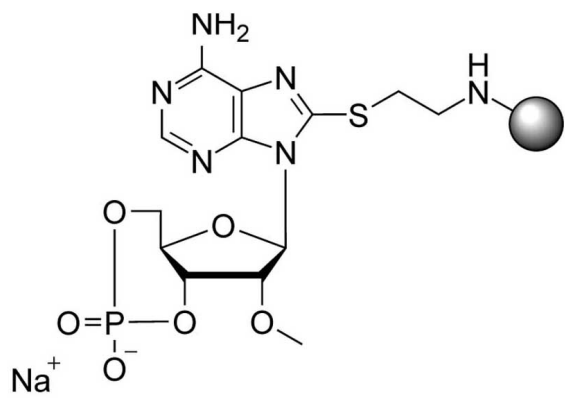
Ordering Information

[Order Online »](#)

BLG-A195-06	0.6ml
BLG-A195-25	2.5ml
BLG-A195-60	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-(2-Aminoethylthio)-2'-O-methyladenosine-3',5'-cyclic monophosphate, 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