

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

8-AET-2'-O-Me-cAMP has only poor affinity towards protein kinases A and G and is thus suitable as a ligand in affinity chromatography of cAMP binding proteins that do not require an intact 2'-OH group, such as the exchange protein activated by cyclic AMP (Epac) and certain phosphodiesterases.

*BLG-A142-25 (5 x 5 µmol pack size) is not sold in the U.S. or Canada.
Please [contact us](#) for available options.*

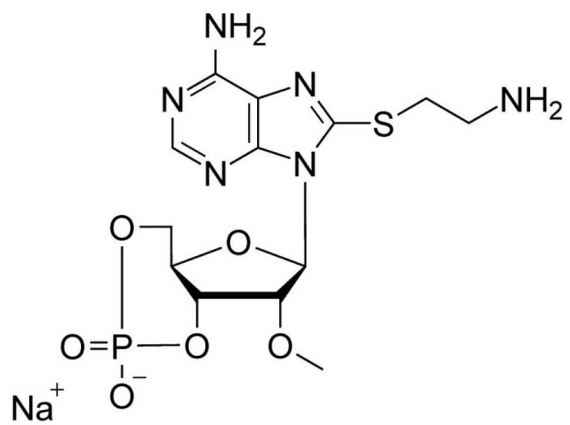
Ordering Information

[Order Online »](#)

BLG-A142-05	5µmol
BLG-A142-25	5x5µmol

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-(2-Aminoethylthio)-2'-O-methyladenosine-3',5'-cyclic monophosphate . sodium salt

CAS 634207-89-9

Couple Type Modified nucleotides

Formula $C_{13}H_{18}N_6O_6PS \cdot Na$

MW 440.4

Purity > 98% HPLC



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