

8-MBT-cAMP

Functionalized cAMP analogue, suitable for immobilization as affinity ligand (e.g. for purification of phosphodiesterases) or for coupling of various labelling structures including fluorophores.

*BLG-M080-25 (5 x 5 μ mol pack size) is not sold in the U.S. or Canada.
Please [contact us](#) for available options.*

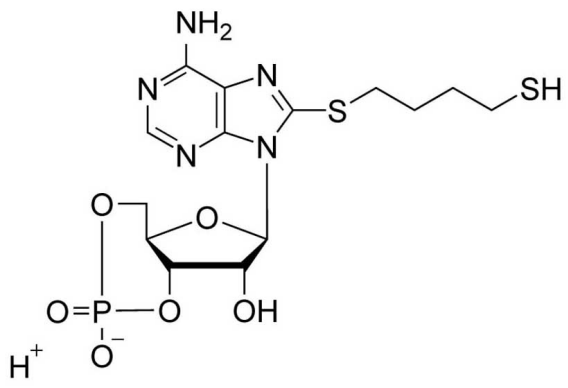
Ordering Information

[Order Online »](#)

BLG-M080-05	5 μ mol
BLG-M080-25	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-(4-Mercaptobutylthio)adenosine-3',5'-cyclic monophosphate . triethyl ammonium salt

CAS 1033893-75-2

Couple Type Modified nucleotides

Formula $C_{14}H_{20}N_5O_6PS_2$ (free acid)

MW 449.4

Purity > 98% HPLC

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