

8-Bromoadenosine 3',5'-cyclic Monophosphothioate, Rp-Isomer . sodium salt

inhibitor of cAMP-dependent protein kinase (PKA), preferentially binding to type I PKA. Lipophilic analog of Rp-cAMPS (Prod. No. BML-CN135, and Prod. No. ALX-480-085). Resistant to mammalian cyclic nucleotide-dependent phosphodiesterases. Significantly more lipophilic and membrane permeant as compared to Rp-cAMPS or 8-Br-cAMP (Prod. No. BML-CN115).

PKA inhibitor

Citations: 3

[View Online »](#)

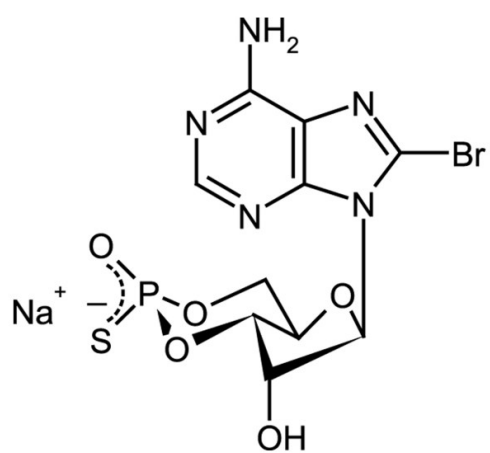
Ordering Information

[Order Online »](#)

ALX-480-024-MC05	0.5mg
ALX-480-024-M001	1mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Solutions should be stored in the refrigerator and should be lyophilized and frozen for longer storage periods.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Rp-8-Br-cAMPS . Na, 8-Br-cAMPS . Na, Rp-isomer
Appearance	White to off-white powder.
CAS	129735-00-8
Couple Target	PKA
Couple Type	Inhibitor
Formula	$C_{10}H_{10}BrN_5O_5PS \cdot Na$
Identity	Identity determined by MS and UV.
MW	423.2 . 23.0
Purity	≥99% (HPLC)
Solubility	250 mM soluble in water, aqueous solutions, DMSO or dimethyl formamide.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)