

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

Metabolically resistant activator of cAMP-dependent protein kinase (PKA) (cAMP agonist). No metabolic side effects as observed with dibutyryl-cAMP (N⁶,O^{2'}-Dibutyryl-adenosine 3',5'-cyclic monophosphate, Prod. No. BML-CN125) or 8-Br-cAMP (8-Bromoadenosine-3',5'-cyclic monophosphate, Prod. No. BML-CN115). Significantly more lipophilic and membrane permeant as compared to Sp-cAMPS (Adenosine 3',5'-cyclic monophosphothioate, Sp isomer, Prod. No. BML-CN136).

Citations: 1

[View Online »](#)

Ordering Information

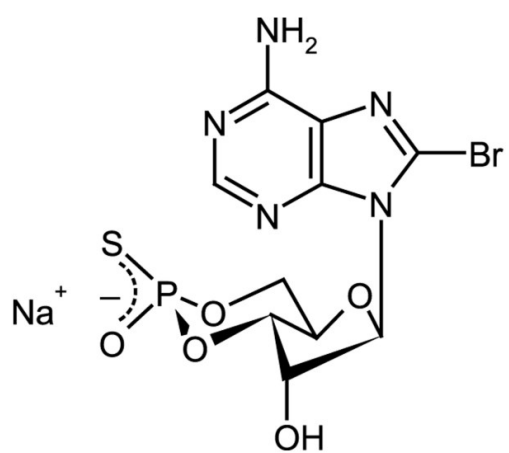
[Order Online »](#)

ALX-480-025-M001

1mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Sp-8-Br-cAMPS . Na, 8-Br-cAMPS . Na, Sp-isomer

Appearance White to off-white powder.

CAS 127634-20-2

Couple Target PKA

Couple Type Activator

Formula $C_{10}H_{10}BrN_5O_5PS \cdot Na$

MW 423.2 . 23.0

Purity ≥98% (HPLC)

Solubility Soluble in water or aqueous buffer.

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)