

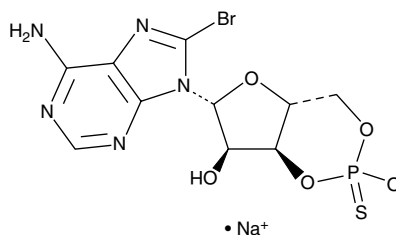
Product Information



Sp-8-bromo-Cyclic AMPS (sodium salt)

Item No. 16002

CAS Registry No.: 1573115-90-8
Formal Name: (4aR,6R,7R,7aS)-6-(6-amino-8-bromo-9H-purin-9-yl)-7-hydroxytetrahydro-4H-furo[3,2-d][1,3,2]dioxaphosphinin-2-olate 2-sulfide, monosodium salt
Synonyms: 8-Bromoadenosine 3',5'-cyclic Monophosphothioate SP-Isomer, Sp-8-bromo-cAMPS
MF: C₁₀H₁₀BrN₅O₅PS • Na
FW: 446.2
Purity: ≥98%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid
UV/Vis.: λ_{max}: 212, 264 nm



Laboratory Procedures

For long term storage, we suggest that Sp-8-bromo-cyclic AMPS (Sp-8-bromo-cAMPS) (sodium salt) be stored as supplied at -20°C. It should be stable for at least two years.

Sp-8-bromo-cAMPS (sodium salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the Sp-8-bromo-cAMPS (sodium salt) in the solvent of choice. Sp-8-bromo-cAMPS (sodium salt) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of Sp-8-bromo-cAMPS (sodium salt) in these solvents is approximately 0.5, 25, and 30 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of Sp-8-bromo-cAMPS (sodium salt) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of Sp-8-bromo-cAMPS (sodium salt) in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Sp-8-bromo-cAMPS is a cell-permeable, cAMP analog that combines an exocyclic sulfur substitution in the axial position of the cyclophosphate ring with a bromine substitution in the adenine base of cAMP.¹⁻² This configuration pools the structural features of two established cyclic-AMP-dependent protein kinase (PKA) activators, 8-bromo-cAMP (Item No. 14431) and Sp-cAMPS (Item No. 14983). Sp-8-bromo-cAMPS is a PKA agonist (EC₅₀ = 1.5 μM) with improved lipophilicity and is not readily degraded by cyclic nucleotide phosphodiesterases.³⁻⁴

References

1. Yokozaki, H., Tortora, G., Pepe, S., *et al.* Unhydrolyzable analogues of adenosine 3':5'-monophosphate demonstrating growth inhibition and differentiation in human cancer cells. *Cancer Res.* **52(9)**, 2504-2508 (1992).
2. Dostmann, W.R., Taylor, S.S., Genieser, H.G., *et al.* Probing the cyclic nucleotide binding sites of cAMP-dependent protein kinases I and II with analogs of adenosine 3',5'-cyclic phosphorothioates. *J. Biol. Chem.* **265(18)**, 10484-10491 (1990).
3. Schwede, F., Maronde, F., Genieser, H., *et al.* Cyclic nucleotide analogs as biochemical tools and prospective drugs. *Pharmacol. Ther.* **87(2)**, 199-226 (2000).
4. Schaap, P., van Ments-Cohen, M., Soede, R.D., *et al.* Cell-permeable non-hydrolyzable cAMP derivatives as tools for analysis of signaling pathways controlling gene regulation in *Dictyostelium*. *J. Biol. Chem.* **268(9)**, 6323-6331 (1993).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/16002

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes no warranty or guarantee of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman warrants only to the original customer that the material will meet our specifications at the time of delivery.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have any obligation or liability, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's exclusive remedy and Cayman's sole liability hereunder shall be limited to a refund of the purchase price, or at Cayman's option, the replacement, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our Warranty and Limitation of Remedy located on our website and in our catalog.

Copyright Cayman Chemical Company, 07/03/2014

Cayman Chemical

Mailing address
1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone
(800) 364-9897
(734) 971-3335

Fax
(734) 971-3640

E-Mail
custserv@caymanchem.com

Web
www.caymanchem.com