

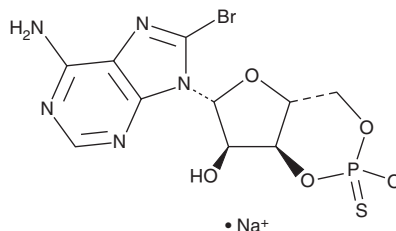
# PRODUCT INFORMATION



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

Item No. 16002

**CAS Registry No.:** 1573115-90-8  
**Formal Name:** cyclic 3',5'-[hydrogen [P(S)]-phosphorothioate]8-bromo-adenosine, monosodium salt  
**Synonyms:** 8-Bromoadenosine 3',5'-cyclic Monophosphothioate SP-Isomer, Sp-8-bromo-cAMPS  
**MF:** C<sub>10</sub>H<sub>10</sub>BrN<sub>5</sub>O<sub>5</sub>PS • Na  
**FW:** 446.2  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 212, 264 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

Sp-8-bromo-cyclic AMPS (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, which should be purged with an inert gas. Sp-8-bromo-cAMPS (sodium salt) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. 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.

### Description

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.<sup>1,2</sup> 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 (EC<sub>50</sub> = 1.5 μM) with improved lipophilicity and is not readily degraded by cyclic nucleotide phosphodiesterases.<sup>3,4</sup>

### References

1. Yokozaki, H., Tortora, G., Pepe, S., et al. *Cancer Res.* **52(9)**, 2504-2508 (1992).
2. Dostmann, W.R., Taylor, S.S., Genieser, H.G., et al. *J. Biol. Chem.* **265(18)**, 10484-10491 (1990).
3. Schwede, F., Maronde, F., Genieser, H., et al. *Pharmacol. Ther.* **87(2)**, 199-226 (2000).
4. Schaap, P., van Ments-Cohen, M., Soede, R.D., et al. *J. Biol. Chem.* **268(9)**, 6323-6331 (1993).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. 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

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/10/2022

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897

[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM