

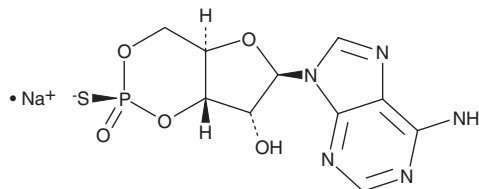
PRODUCT INFORMATION



Sp-Cyclic AMPS (sodium salt)

Item No. 14983

CAS Registry No.: 142439-95-0
Formal Name: cyclic 3',5'-[hydrogen (S)-phosphorothioate] adenosine, monosodium salt
Synonym: Sp-cAMPS
MF: C₁₀H₁₁N₅O₅PS • Na
FW: 367.3
Purity: ≥98%
UV/Vis.: λ_{max}: 259 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-cAMPS (sodium salt) is supplied as a crystalline solid. Sp-cAMPS (sodium salt) is soluble in water at a concentration of approximately 25 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Several cyclic nucleotide phosphodiesterases (PDE2, -5, -6, -10, and -11) contain a tandem of GAF domains in their N-terminal regulatory region that are thought to mediate cAMP- or cGMP-induced stimulation of adenylyl and guanylyl cyclases.¹ Sp-cAMPS is a derivative of cAMP that binds the PDE10 GAF domain (EC₅₀ = 40 μM).¹ It has been used to study GAF-mediated stimulation of PDE10.¹

Reference

1. Jäger, R., Russwurm, C., Schwede, F., *et al.* Activation of PDE10 and PDE11 phosphodiesterases. *J. Biol. Chem.* **287**(2), 1210-1219 (2012).

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, 11/23/2022

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