

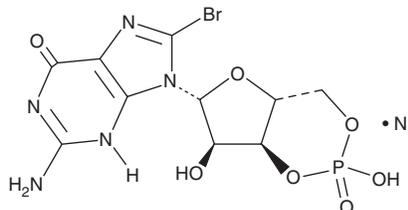
PRODUCT INFORMATION



8-bromo Cyclic GMP (sodium salt)

Item No. 15992

CAS Registry No.: 51116-01-9
Formal Name: 8-bromo-guanosine cyclic 3',5'-(hydrogen phosphate), monosodium salt
Synonyms: 8-bromo cGMP, 8-bromo Guanosine 3',5'-cyclic monophosphate
MF: C₁₀H₁₁BrN₅O₇P • Na
FW: 447.1
Purity: ≥98%
UV/Vis.: λ_{max}: 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

8-bromo cGMP (sodium salt) is supplied as a crystalline solid. Aqueous solutions of 8-bromo cGMP (sodium salt) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of 8-bromo cGMP (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

8-bromo cGMP is a cell-permeable analog of cGMP that exhibits greater resistance to hydrolysis by phosphodiesterases than its parent compound. It preferentially activates cGMP-dependent protein kinase.¹ At 1 μM-0.1 mM, 8-bromo cGMP can inhibit acetylcholine-induced increases in intracellular calcium concentrations.²

References

1. Rashatwar, S.S., Cornwell, T.L., and Lincoln, T.M. Effects of 8-bromo-cGMP on Ca²⁺ levels in vascular smooth muscle cells: Possible regulation of Ca²⁺-ATPase by cGMP-dependent protein kinase. *Proc. Natl. Acad. Sci. USA* **84(16)**, 5685-5689 (1987).
2. Choi, J. and Farley, J.M. Effects of 8-bromo-cyclic GMP on membrane potential of single swine tracheal smooth muscle cells. *J. Pharmacol. Exp. Ther.* **285(2)**, 588-594 (1998).

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, 07/05/2023

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