

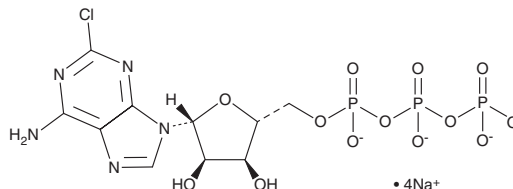
PRODUCT INFORMATION



2-Chloroadenosine 5-triphosphate (sodium salt)

Item No. 34978

CAS Registry No.: 301334-89-4
Formal Name: 2-chloro-adenosine
5'-(tetrahydrogen triphosphate),
tetrasodium salt
Synonyms: 2-chloro ATP
MF: C₁₀H₁₁ClN₅O₁₃P₃ • 4Na
FW: 629.6
Purity: ≥95%
Supplied as: A solution in water
Storage: -80°C
Stability: ≥2 years



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

Description

2-Chloroadenosine 5-triphosphate (2-chloro ATP) is an adenine nucleotide and analog of ATP (Item No. 14498).^{1,2} It is an antagonist of the purinergic P2Y₁ receptor, inhibiting intracellular calcium mobilization induced by ADP (Item No. 21121) in Jurkat cells expressing the human receptor (K_i = 2.3 μM).³ 2-chloro ATP is an agonist of purinergic P2X receptors, inducing inward currents in HEK293 cells expressing the human bladder smooth muscle or rat PC12 forms of the receptor (EC₅₀s = 0.5 and 2.5 μM, respectively).⁴ It induces relaxation of precontracted isolated guinea pig taenia caeci strips in a concentration-dependent manner.¹ It has been used to study the substrate specificity of cyclic nucleotide-dependent protein kinases, such as protein kinase A (PKA) and PKG.²

References

1. Satchell, G. and Maguire, M.H. Inhibitory effects of adenine nucleotide analogs on the isolated guinea-pig taenia coli. *J. Pharmacol. Exp. Ther.* **195**(3), 540-548 (1975).
2. Flockhart, D.A., Freist, W., Hoppe, J.O., et al. ATP analog specificity of cAMP-dependent protein kinase, cGMP-dependent protein kinase, and phosphorylase kinase. *Eur. J. Biochem.* **140**(2), 289-295 (1984).
3. Hechler, B., Vigne, P., Léon, C., et al. ATP derivatives are antagonists of the P2Y1 receptor: Similarities to the platelet ADP receptor. *Mol. Pharmacol.* **53**(4), 727-733 (1997).
4. Evans, R.J., Lewis, C., Buell, G., et al. Pharmacological characterization of heterologously expressed ATP-gated cation channels (P2x purinoceptors). *Mol. Pharmacol.* **48**(2), 178-183 (1995).

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, 03/30/2023

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