

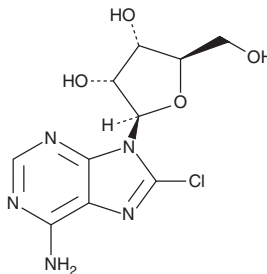
# PRODUCT INFORMATION



## 8-Chloroadenosine

Item No. 35766

CAS Registry No.: 34408-14-5  
Formal Name: 8-chloro-adenosine  
MF:  $C_{10}H_{12}ClN_5O_4$   
FW: 301.7  
Purity:  $\geq 98\%$   
UV/Vis.:  $\lambda_{max}$ : 262 nm  
Supplied as: A solid  
Storage:  $-20^{\circ}C$   
Stability:  $\geq 4$  years



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

### Laboratory Procedures

8-Chloroadenosine is supplied as a solid. A stock solution may be made by dissolving the 8-chloroadenosine in the solvent of choice, which should be purged with an inert gas. 8-Chloroadenosine is slightly soluble in DMSO.

### Description

8-Chloroadenosine is a nucleoside analog and an active metabolite of the anticancer agent 8-chloro cAMP.<sup>1</sup> It is formed from 8-chloro cAMP by phosphodiesterases (PDEs) and upon entry into cells, 8-chloroadenosine is phosphorylated by adenosine kinase to produce 8-chloro-ATP. 8-Chloroadenosine (10  $\mu$ M) inhibits RNA, but not DNA, synthesis *in vitro*. It is cytotoxic to patient-derived multiple myeloma cells in a concentration-dependent manner. 8-Chloroadenosine (50 mg/kg) reduces tumor growth by 50% in an HCT116 colorectal cancer mouse xenograft model.<sup>2</sup>

### References

1. Gandhi, V., Ayres, M., Halgren, R.G., *et al.* 8-chloro-cAMP and 8-chloro-adenosine act by the same mechanism in multiple myeloma cells. *Cancer Res.* **61(14)**, 5474-5479 (2001).
2. Carlson, C.C., Chinery, R., Burnham, L.L., *et al.* 8-Cl-adenosine-induced inhibition of colorectal cancer growth *in vitro* and *in vivo*. *Neoplasia* **2(5)**, 441-448 (2000).

#### 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, 12/13/2022

#### 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