

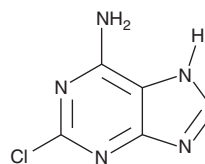
# PRODUCT INFORMATION



## 2-Chloroadenine

Item No. 30474

**CAS Registry No.:** 1839-18-5  
**Formal Name:** 2-chloro-9H-purin-6-amine  
**Synonyms:** 6-amino-2-Chloropurine, 2-chloro-6-Aminopurine, NSC 7362  
**MF:** C<sub>5</sub>H<sub>4</sub>ClN<sub>5</sub>  
**FW:** 169.6  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 265 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥2 years



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

### Laboratory Procedures

2-Chloroadenine is supplied as a crystalline solid. A stock solution may be made by dissolving the 2-chloroadenine in the solvent of choice, which should be purged with an inert gas. 2-Chloroadenine is slightly soluble in ethanol.

### Description

2-Chloroadenine is a heterocyclic building block that has been used in the synthesis of adenosine antimetabolites with antineoplastic activity.<sup>1</sup> It is also the major catabolite of 2-chloro-2'-deoxyadenosine.<sup>2</sup>

### References

1. Bauta, W.E., Schulmeier, B.E., Burke, B.J., *et al.* A new process for antineoplastic agent clofarabine. *Org. Proc. Res. Dev.* **8(6)**, 889-896 (2004).
2. Bontemps, F., Delacauw, A., Cardoen, S., *et al.* Metabolism and cytotoxic effects of 2-chloroadenine, the major catabolite of 2-chloro-2'-deoxyadenosine. *Biochem. Pharmacol.* **59(10)**, 1237-1243 (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, 05/04/2020

#### 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](http://WWW.CAYMANCHEM.COM)