

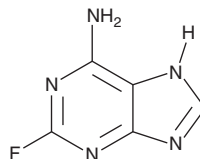
# PRODUCT INFORMATION



## 2-Fluoroadenine

Item No. 30421

CAS Registry No.: 700-49-2  
Formal Name: 2-fluoro-9H-purin-6-amine  
Synonyms: F-Ade, NSC 27364  
MF:  $C_5H_4FN_5$   
FW: 153.1  
Purity:  $\geq 98\%$   
UV/Vis.:  $\lambda_{\max}$ : 262 nm  
Supplied as: A crystalline solid  
Storage:  $-20^{\circ}\text{C}$   
Stability:  $\geq 4$  years



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

### Laboratory Procedures

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

### Description

2-Fluoroadenine is a heterocyclic building block that has been used in the synthesis of heat shock protein 90 (Hsp90) inhibitors that have anticancer activity *in vitro*.<sup>1</sup> It is also cytotoxic to CEM human leukemia cells with an  $IC_{50}$  value of  $0.15\ \mu\text{M}$ .<sup>2</sup>

### References

1. Llauger, L., He, H., Kim, J., *et al.* Evaluation of 8-arylsulfanyl, 8-arylsulfoxyl, and 8-arylsulfonyl adenine derivatives as inhibitors of the heat shock protein 90. *J. Med. Chem.* **48**(8), 2892-2905 (2005).
2. Parker, W.B., Allan, P.W., Shaddix, S.C., *et al.* Metabolism and metabolic actions of 6-methylpurine and 2-fluoroadenine in human cells. *Biochem. Pharmacol.* **55**(10), 1673-1681 (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, 11/22/2022

#### 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