

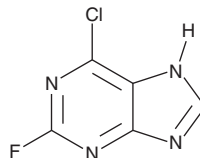
PRODUCT INFORMATION



6-Chloro-2-fluoropurine

Item No. 30481

CAS Registry No.: 1651-29-2
Synonym: NSC 37363
MF: $C_5H_2ClFN_4$
FW: 172.5
Purity: $\geq 98\%$
UV/Vis.: λ_{\max} : 270 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

6-Chloro-2-fluoropurine is supplied as a crystalline solid. A stock solution may be made by dissolving the 6-chloro-2-fluoropurine in the solvent of choice, which should be purged with an inert gas. 6-Chloro-2-fluoropurine is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of 6-chloro-2-fluoropurine in these solvents is approximately 30 mg/ml.

Description

6-Chloro-2-fluoropurine is a heterocyclic building block.^{1,2} It has been used in the synthesis of purine nucleosides that inhibit cyclin-dependent kinases (CDKs) *in vitro*.¹ 6-Chloro-2-fluoropurine has also been used in the synthesis of purine nucleosides that are active against HIV-1 and hepatitis B virus (HBV) *in vitro*.²

References

1. Wilson, S.C., Atrash, B., Barlow, C., *et al.* Design, synthesis and biological evaluation of 6-pyridylmethylaminopurines as CDK inhibitors. *Bioorg. Med. Chem.* **19(22)**, 6949-6965 (2011).
2. Lee, K., Choi, Y., Gullen, E., *et al.* Synthesis and anti-HIV and anti-HBV activities of 2'-fluoro-2',3'-unsaturated L-nucleosides. *J. Med. Chem.* **42(7)**, 1320-1328 (1999).

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, 11/03/2022

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