

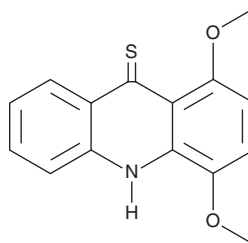
PRODUCT INFORMATION



NSC 625987

Item No. 32843

CAS Registry No.: 141992-47-4
Formal Name: 1,4-dimethoxy-9(10H)-acridinethione
Synonym: Cdk4 Inhibitor II
MF: C₁₅H₁₃NO₂S
FW: 271.3
Purity: ≥98%
Supplied as: A 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

NSC 625987 is supplied as a solid. A stock solution may be made by dissolving the NSC 625987 in the solvent of choice, which should be purged with an inert gas. NSC 625987 is soluble in ethanol and DMSO.

Description

NSC 625987 is an inhibitor of the Cdk4/cyclin D1 complex ($IC_{50} = 0.2 \mu M$).^{1,2} It is selective for the Cdk4/cyclin D1 complex over Cdk2/cyclin A, Cdk2/cyclin E, and Cdk1/cyclin A complexes (IC_{50} s = >100 μM for all).¹

References

1. Kubo, A., Nakagawa, K., Varma, R.K., *et al.* The p16 status of tumor cell lines identifies small molecule inhibitors specific for cyclin-dependent kinase 4. *Clin. Cancer Res.* **5(12)**, 4279-4286 (1999).
2. Jorda, R., Hendrychová, D., Voller, J., *et al.* How selective are pharmacological inhibitors of cell-cycle-regulating cyclin-dependent kinases? *J. Med. Chem.* **61(20)**, 9105-9120 (2018).

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, 01/03/2023

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