

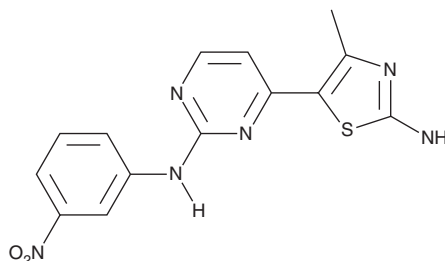
PRODUCT INFORMATION



CK7

Item No. 39153

CAS Registry No.: 507487-89-0
Formal Name: 4-(2-amino-4-methyl-5-thiazolyl)-N-(3-nitrophenyl)-2-pyrimidinamine
Synonyms: Cdk2/9 Inhibitor, Cyclin-dependent kinase 2/9 Inhibitor
MF: C₁₄H₁₂N₆O₂S
FW: 328.4
Purity: ≥95%
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

CK7 is supplied as a solid. A stock solution may be made by dissolving the CK7 in the solvent of choice, which should be purged with an inert gas. CK7 is slightly soluble in methanol and DMSO.

Description

CK7 is an inhibitor of cyclin-dependent kinase 2 (Cdk2) and Cdk9 (K_i s = 2 and 4 nM, respectively).¹ It is selective for Cdk2 and Cdk9 over Cdk1, Cdk4, and Cdk7 (K_i s = 80, 53, and 70 nM, respectively), as well as 10 additional kinases (K_i s = >1,000 nM for all). CK7 is cytotoxic to A2780 ovarian and MES-SA uterine cancer cells (IC_{50} s = 41 and 140 nM, respectively).²

References

1. Wang, S., Meades, C., Wood, G., *et al.* 2-Anilino-4-(thiazol-5-yl)pyrimidine CDK inhibitors: Synthesis, SAR analysis, X-ray crystallography, and biological activity. *J. Med. Chem.* **47**(7), 1662-1675 (2004).
2. Wang, S., Griffiths, G., Midgley, C.A., *et al.* Discovery and characterization of 2-anilino-4-(thiazol-5-yl)pyrimidine transcriptional CDK inhibitors as anticancer agents. *Chem. Biol.* **17**(10), 1111-1121 (2010).

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, 12/12/2023

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