

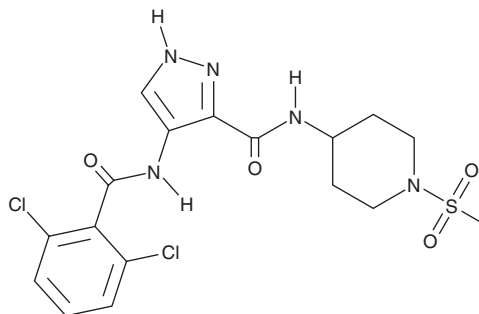
PRODUCT INFORMATION



NVP-LCQ195

Item No. 42728

CAS Registry No.: 902156-99-4
Formal Name: 4-[(2,6-dichlorobenzoyl)amino]-N-[1-(methylsulfonyl)-4-piperidiny]-1H-pyrazole-3-carboxamide
Synonyms: AT-9311, LCQ-195
MF: C₁₇H₁₉Cl₂N₅O₄S
FW: 460.3
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

NVP-LCQ195 is supplied as a solid. A stock solution may be made by dissolving the NVP-LCQ195 in the solvent of choice, which should be purged with an inert gas. NVP-LCQ195 is slightly soluble (0.1-1 mg/ml) in ethanol.

Description

NVP-LCQ195 is an inhibitor of cyclin-dependent kinases (CDKs).¹ It selectively inhibits Cdk5/p25, Cdk5/p35, Cdk1/cyclin B, Cdk2/cyclin A, and Cdk2/cyclin E complexes (IC₅₀s = 1, 1, 2, 2, and 5 nM, respectively) over Cdk9/cyclin T1, Cdk3/cyclin E, Cdk6/cyclin D3, and Cdk7/cyclin H complexes (IC₅₀s = 15, 42, 187, and 3,564 nM, respectively). NVP-LCQ195 decreases viability in a panel of 25 multiple myeloma cells (EC₅₀s = 0.761-2.81 μM). It induces cell cycle arrest at the G₀ phase and apoptosis in MM.1S multiple myeloma cells when used at a concentration of 2 μM.

Reference

1. McMillin, D.W., Delmore, J., Negri, J., *et al.* Molecular and cellular effects of multi-targeted cyclin-dependent kinase inhibition in myeloma: Biological and clinical implications. *Br. J. Haematol.* **152**(4), 420-432 (2011).

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, 02/07/2025

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