

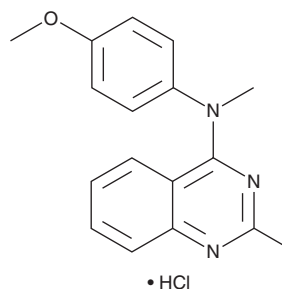
PRODUCT INFORMATION



MPC-6827 (hydrochloride)

Item No. 32872

CAS Registry No.: 917369-31-4
Formal Name: N-(4-methoxyphenyl)-N,2-dimethyl-4-quinazolinamine, monohydrochloride
Synonym: Verubulin
MF: C₁₇H₁₇N₃O • HCl
FW: 315.8
Purity: ≥98%
UV/Vis.: λ_{max}: 223, 334 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

MPC-6827 (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the MPC-6827 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. MPC-6827 (hydrochloride) is soluble in the organic solvent DMSO. It is also soluble in water. The solubility of MPC-6827 (hydrochloride) in DMSO and water is approximately 100 mM. We do not recommend storing the aqueous solution for more than one day.

Description

MPC-6827 is an inhibitor of tubulin polymerization with anticancer activity.^{1,2} It inhibits tubulin polymerization in a cell-free assay when used at concentrations of 0.5, 5, and 50 μM, as well as in A549 cells when used at a concentration of 10 nM.¹ MPC-6827 induces apoptosis in T47D, HCT116, and SNU-398 cancer cells (EC₅₀s = 2, 3, and 2 nM, respectively).² *In vivo*, MPC-6827 (2.5-10 mg/kg) reduces tumor growth in MX-1, OVCAR-3, MIA PaCa-2, MCF-7, HT-29, and MDA-MB-435 mouse xenograft models.¹

References

1. Kasibhatla, S., Baichwal, V., Cai, S.X., *et al.* MPC-6827: A small-molecule inhibitor of microtubule formation that is not a substrate for multidrug resistance pumps. *Cancer Res.* **67(12)**, 5865-5871 (2007).
2. Sirisoma, N., Pervin, A., Zhang, H., *et al.* Discovery of N-(4-methoxyphenyl)-N,2-dimethylquinazolin-4-amine, a potent apoptosis inducer and efficacious anticancer agent with high blood brain barrier penetration. *J. Med. Chem.* **52(8)**, 2341-2351 (2009).

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, 11/10/2022

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