

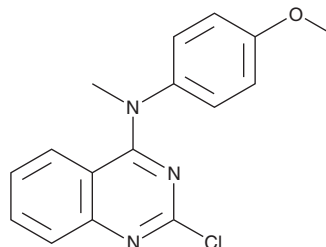
PRODUCT INFORMATION



MPI-0441138

Item No. 13098

CAS Registry No.: 827030-33-1
Formal Name: 2-chloro-N-(4-methoxyphenyl)-N-methyl-4-quinazolinamine
Synonym: EP128265
MF: C₁₆H₁₄ClN₃O
FW: 299.8
Purity: ≥98%
UV/Vis.: λ_{max}: 215, 383, 336 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

MPI-0441138 is supplied as a crystalline solid. A stock solution may be made by dissolving the MPI-0441138 in the solvent of choice, which should be purged with an inert gas. MPI-0441138 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of MPI-0441138 in ethanol is approximately 5 mg/ml and approximately 10 mg/ml in DMSO and DMF.

MPI-0441138 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, MPI-0441138 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. MPI-0441138 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

MPI-0441138 is a potent inducer of apoptosis and growth inhibition (EC₅₀ = 2 nM for both processes, based on caspase-3 activation and total cellular ATP, respectively) in T47D and HCT116 cells.^{1,2} MPI-0441138 has excellent blood brain barrier penetration and is effective in mouse xenograft cancer models.^{1,2}

References

1. Sirisoma, N., Kasibhatla, S., Pervin, A., *et al.* Discovery of 2-chloro-N-(4-methoxyphenyl)-N-methylquinazolin-4-amine (EP128265, MPI-0441138) as a potent inducer of apoptosis with high *in vivo* activity. *J. Med. Chem.* **51**, 4771-4779 (2008).
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, 11/22/2022

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