

PRODUCT INFORMATION



M606

Item No. 45735

CAS Registry No.: 298230-09-8

Formal Name: 5-(methoxymethyl)-7-[(4-methyl-1-piperidinyl)methyl]-8-quinolinol

MF: C₁₈H₂₄N₂O₂

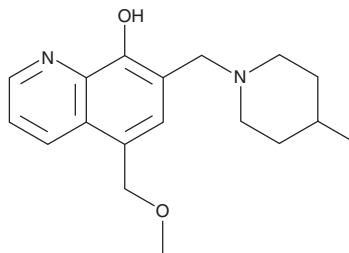
FW: 300.4

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

M606 is supplied as a solid. A stock solution may be made by dissolving the M606 in the solvent of choice, which should be purged with an inert gas. M606 is soluble (≥10 mg/ml) in ethanol and sparingly soluble (1-10 mg/ml) in DMSO.

Description

M606 is an iron chelator and N-Myc inhibitor.¹ It binds iron at a 3:1 ratio. M606 inhibits MYCN transactivation in a reporter assay using SHR6-17 neuroblastoma cells in a concentration-dependent manner. It induces cytotoxicity, decreases N-Myc levels, and increases HIF-1α levels in BE(2)-C and LA-N-1 neuroblastoma cells, effects that can be reversed by ferric iron. M606 (25 mg/kg) increases survival in the *TH-MYCN* transgenic mouse model of neuroblastoma. It also inhibits *P. aeruginosa* biofilm formation (EC₅₀ = 19.21 μM).²

References

1. Pandher, R., Xue, C., Gamble, L.D., *et al.* The cell-permeable iron chelator M606 inhibits MYCN-driven neuroblastoma via an E2F3-mediated response. *Proc. Nat. Acad. Sci. USA* **122(23)**, e2420011122 (2025).
2. Junker, L.M. and Clardy, J. High-throughput screens for small-molecule inhibitors of *Pseudomonas aeruginosa* biofilm development. *Antimicrob. Agents Chemother.* **51(10)**, 3582-3590 (2007).

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, 07/09/2026

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