

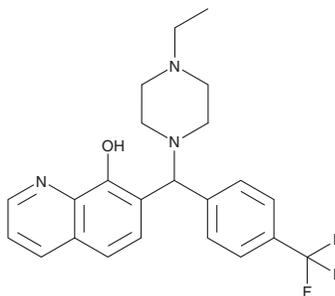
PRODUCT INFORMATION



ML-311

Item No. 29445

CAS Registry No.: 315698-17-0
Formal Name: 7-[(4-ethyl-1-piperazinyl)
[4-(trifluoromethyl)phenyl]
methyl]-8-quinolinol
MF: C₂₃H₂₄F₃N₃O
FW: 415.5
Purity: ≥98%
UV/Vis.: λ_{max}: 250 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

ML-311 is supplied as a crystalline solid. A stock solution may be made by dissolving the ML-311 in the solvent of choice, which should be purged with an inert gas. ML-311 is soluble in the organic solvent DMSO.

Description

ML-311 is an inhibitor of myeloid cell leukemia-1 (Mcl-1; IC₅₀ = 0.31 μM in a fluorescence polarization assay).¹ It is selective for Mcl-1 over Bcl-xL (IC₅₀ = >40 μM). ML-311 decreases the proliferation of 13 cancer cell lines (EC₅₀s = 0.3-25 μM).

Reference

1. Richard, D.J., Lena, R., Bannister, T., *et al.* Hydroxyquinoline-derived compounds and analoguing of selective Mcl-1 inhibitors using a functional biomarker. *Bioorg. Med. Chem.* **21(21)**, 6642-6649 (2013).

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, 11/04/2022

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