

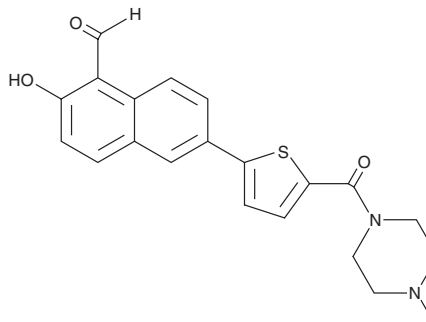
PRODUCT INFORMATION



MKC-3946

Item No. 19152

CAS Registry No.: 1093119-54-0
Formal Name: 2-hydroxy-6-[5-[(4-methyl-1-piperazinyl)carbonyl]-2-thienyl]-1-naphthalenecarboxaldehyde
MF: C₂₁H₂₀N₂O₃S
FW: 380.5
Purity: ≥98%
UV/Vis.: λ_{max}: 226, 316 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

MKC-3946 is supplied as a crystalline solid. A stock solution may be made by dissolving the MKC-3946 in the solvent of choice, which should be purged with an inert gas. MKC-3946 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of MKC-3946 in these solvents is approximately 14 and 20 mg/ml, respectively.

MKC-3946 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, MKC-3946 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. MKC-3946 has a solubility of approximately 0.14 mg/ml in a 1:6 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

MKC-3946 is an inhibitor of inositol-requiring enzyme 1α (IRE1α).¹ It inhibits basal XBP1 splicing, a marker of IRE1α activity, in RPMI 8226 multiple myeloma cells in a concentration-dependent manner. MKC-3946 also inhibits XBP1 splicing induced by endoplasmic reticulum (ER) stress and reduces expression of the XBP1 target genes SEC61A1, p58IPK, and ERdj4 in RPMI 8226 cells. *In vitro*, it is cytotoxic to multiple myeloma cell lines but has no effect on normal mononuclear cells. *In vivo*, MKC-3946 (100 mg/kg) reduces tumor growth and XBP1 splicing in an RPMI 8226 mouse xenograft model.

Reference

1. Mimura, N., Fulciniti, M., Gorgun, G., *et al.* Blockade of XBP1 splicing by inhibition of IRE1α is a promising therapeutic option in multiple myeloma. *Blood* **119(24)**, 5772-5781 (2012).

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, 10/18/2022

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