

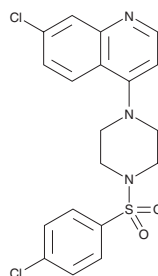
PRODUCT INFORMATION



KM 11060

Item No. 23481

CAS Registry No.: 774549-97-2
Formal Name: 7-chloro-4-[4-[(4-chlorophenyl)sulfonyl]-1-piperazinyl]-quinoline
MF: C₁₉H₁₇Cl₂N₃O₂S
FW: 422.3
Purity: ≥98%
UV/Vis.: λ_{max}: 201, 227, 321 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

KM 11060 is supplied as a crystalline solid. A stock solution may be made by dissolving the KM 11060 in the solvent of choice, which should be purged with an inert gas. KM 11060 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of KM 11060 in these solvents is approximately 0.5, 2, and 5 mg/ml, respectively.

Description

KM 11060 is a small molecule that corrects the processing of cystic fibrosis transmembrane conductance regulator (CFTR) proteins bearing the F508 deletion (F508del) mutation carried by 90% of cystic fibrosis patients.¹ It increases cell surface expression and restores chloride transport function of F508del-CFTRs in BHK cells at a concentration of 10 μM. KM 11060 restores F508del-CFTR function, as measured by cAMP-stimulated iodide efflux, in CFBE41o⁻ human airway epithelial cells at a concentration of 10 nM. It also restores the secretory response to approximately 51% of wild-type levels *ex vivo* in ileum isolated from F508del-CFTR mice at a concentration of 20 μM.

Reference

1. Robert, R., Carlile, G.W., Pavel, C. *et al.* Structural analog of sildenafil identified as a novel corrector of the F508del-CFTR trafficking defect. *Mol. Pharmacol.* **73(2)**, 478-489 (2008).

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, 11/15/2022

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