

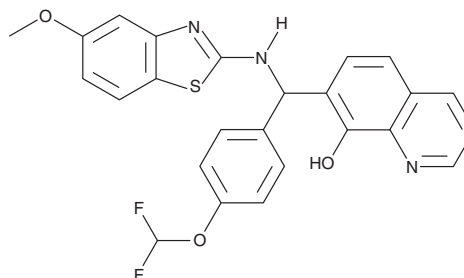
PRODUCT INFORMATION



KIN1408

Item No. 21790

CAS Registry No.: 1903800-11-2
Formal Name: 7-[[4-(difluoromethoxy)phenyl]
[(5-methoxy-2-benzothiazolyl)amino]
methyl]-8-quinolinol
MF: C₂₅H₁₉F₂N₃O₃S
FW: 479.5
Purity: ≥98%
UV/Vis.: λ_{max}: 247, 302, 309 nm
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

KIN1408 is supplied as a solid. A stock solution may be made by dissolving the KIN1408 in the solvent of choice. KIN1408 is soluble in the organic solvent DMSO, which should be purged with an inert gas, at a concentration of approximately 30 mg/ml.

Description

KIN1408 is a small molecule activator of the RIG-1-like receptor (RLR) pathway that has antiviral activity.¹ It induces interferon regulatory factor 3 (IRF3) activation and translocation into the nucleus in a concentration-dependent manner in HEK293 cells without inducing cellular toxicity. KIN1408 induces expression of the innate immune genes MDA5, RIG-1, Mx1, IRF7, and IFIT1 in THP-1 cells stimulated with phorbol 12-myristate 12-acetate (PMA; Item No. 10008014). It exhibits concentration-dependent broad-spectrum antiviral activity against dengue virus 2 as well as the influenza A, Ebola, Nipah, and Lassa viruses.

Reference

1. Pattabhi, S., Wilkins, C.R., Dong, R., *et al.* Targeting innate immunity for antiviral therapy through small molecule agonists of the RLR Pathway. *J. Virol.* **90**(5), 2372-2387 (2015).

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, 12/12/2022

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