

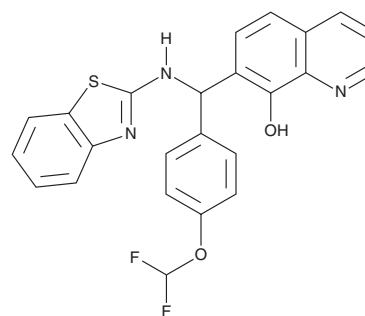
PRODUCT INFORMATION



KIN1400

Item No. 22441

CAS Registry No.: 446826-86-4
Formal Name: 7-[(2-benzothiazolylamino)[4-(difluoromethoxy)phenyl]methyl]-8-quinolinol
MF: C₂₄H₁₇F₂N₃O₂S
FW: 449.5
Purity: ≥98%
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

KIN1400 is supplied as a solid. A stock solution may be made by dissolving the KIN1400 in the solvent of choice, which should be purged with an inert gas. KIN1400 is soluble in the organic solvent DMSO at a concentration of approximately 30 mg/ml.

Description

KIN1400 is a small molecule activator of the RIG-1-like receptor (RLR) pathway that has antiviral activity.¹ It induces expression of the innate immune genes RIG-1, MDA5, IFIT1, IFIT2, IFITM1, OAS3, and Mx1 in THP-1 cells stimulated with phorbol 12-myristate 13-acetate (PMA; Item No. 10008014). KIN1400 (20 µM) reduces West Nile and dengue viral RNA levels in HEK293 and Huh7 cells, respectively. It also inhibits hepatitis C virus (HCV) replication in a concentration-dependent manner in Huh7 cells.

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, 11/16/2022

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