

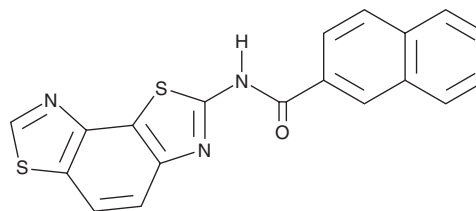
PRODUCT INFORMATION



KIN1148

Item No. 38980

CAS Registry No.: 1428729-56-9
Formal Name: N-benzo[1,2-d:3,4-d']bisthiazol-2-yl-2-naphthalenecarboxamide
MF: C₁₉H₁₁N₃OS₂
FW: 361.4
Purity: ≥98%
UV/Vis.: λ_{max}: 217, 249, 310, 326 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

KIN1148 is supplied as a crystalline solid. A stock solution may be made by dissolving the KIN1148 in the solvent of choice, which should be purged with an inert gas. KIN1148 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of KIN1148 in these solvents is approximately 5 mg/ml. KIN1148 is slightly soluble in ethanol.

Description

KIN1148 is an IFN regulating factor 3 (IRF3) activator and adjuvant.¹ It increases the mRNA expression of the genes encoding IFN-stimulated gene 54 (ISG54) and 2'-5'-oligoadenylate synthase-like (OASL) protein in PH5CH8 cells and increases chemokine (C-X-C motif) ligand 10 (CXCL10) production in phorbol 12-myristate 13-acetate-differentiated THP-1 monocytes when used at a concentration of 10 μM. Immunization with a vaccine in combination with KIN1148 as an adjuvant increases blood IgG, neutralizing, and hemagglutinin antibody titers, decreases particle-forming units (PFUs) in the lungs and lung-draining lymph nodes, and increases the survival percentage of mice infected with a lethal dose of influenza virus.

Reference

1. Probst, P., Grigg, J.B., Wang, M., *et al.* A small-molecule IRF3 agonist functions as an influenza vaccine adjuvant by modulating the antiviral immune response. *Vaccine* **35(15)**, 1964-1971 (2017).

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, 07/14/2023

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