

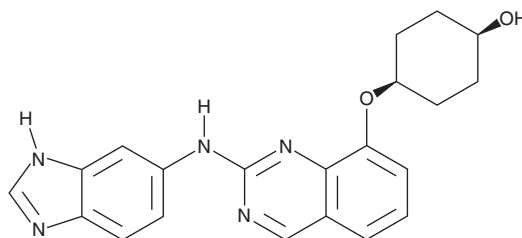
PRODUCT INFORMATION



NCB-0846

Item No. 33268

CAS Registry No.: 1792999-26-8
Formal Name: cis-4-[[2-(1H-benzimidazol-6-ylamino)-8-quinazolinyl]oxy]-cyclohexanol
MF: C₂₁H₂₁N₅O₂
FW: 375.4
Purity: ≥90%
UV/Vis.: λ_{max}: 222, 246, 293, 311 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

NCB-0846 is supplied as a crystalline solid. A stock solution may be made by dissolving the NCB-0846 in the solvent of choice, which should be purged with an inert gas. NCB-0846 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of NCB-0846 in these solvents is approximately 5 and 1 mg/ml, respectively.

Description

NCB-0846 is an inhibitor of TRAF2- and NCK-interacting kinase (TNIK; IC₅₀ = 21 nM).¹ It is selective for TNIK over a panel of 46 kinases but does inhibit FLT3, JAK3, PDGFRα, TrkA, Cdk2/CycA2, and HGK by greater than 80% at 100 nM. NCB-0846 (3 μM) inhibits phosphorylation of transcription factor 4 (TCF4) and autophosphorylation of TNIK in cell-based assays. It inhibits the growth and colony formation of HCT116 colon cancer cells in two-dimensional growth and soft-agar assays, respectively, as well as reduces tumor growth in an HCT116 mouse xenograft model. NCB-0846 (50 and 100 mg/kg) also reduces tumor growth in patient-derived xenograft (PDX) mouse models of colon cancer.

Reference

1. Masuda, M., Uno, Y., Ohbayashi, N., *et al.* TNIK inhibition abrogates colorectal cancer stemness. *Nat. Commun.* **7**, 12586 (2016).

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, 10/17/2022

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