

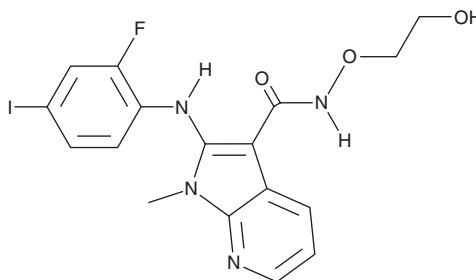
PRODUCT INFORMATION



Nedometinib

Item No. 40465

CAS Registry No.: 2252314-46-6
Formal Name: 2-[(2-fluoro-4-iodophenyl)amino]-N-(2-hydroxyethoxy)-1-methyl-1H-pyrrolo[2,3-b]pyridine-3-carboxamide
Synonym: NFX-179
MF: C₁₇H₁₆FIN₄O₃
FW: 470.2
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

Nedometinib is supplied as a solid. A stock solution may be made by dissolving the nedometinib in the solvent of choice, which should be purged with an inert gas. Nedometinib is soluble in DMSO and methanol.

Description

Nedometinib is a MEK1 inhibitor (IC₅₀ = 135 nM).¹ It is selective for MEK1 over a panel of 44 receptors, ion channels, and enzymes at 10 μM, as well as a panel of an additional 150 kinases. Nedometinib (0.1-10 μM) decreases phosphorylation of ERK in malignant peripheral nerve sheath tumor cells. Topical application of nedometinib (0.1-2.3%) reduces the total number of tumors in a mouse model of cutaneous squamous cell carcinoma induced by UV radiation.

Reference

1. Sarin, K.Y., J., K., B., S., *et al.* Development of a MEK inhibitor, NFX-179, as a chemoprevention agent for squamous cell carcinoma. *Sci. Transl. Med.* **15(717)**, (2023).

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, 03/25/2024

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