

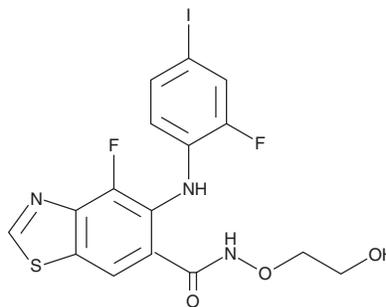
PRODUCT INFORMATION



Tunlametinib

Item No. 45050

CAS Registry No.: 1801756-06-8
Formal Name: 4-fluoro-5-[(2-fluoro-4-iodophenyl)amino]-N-(2-hydroxyethoxy)-6-benzothiazolecarboxamide
Synonym: HL 085
MF: C₁₆H₁₂F₂IN₃O₃S
FW: 491.3
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

Tunlametinib is supplied as a solid. A stock solution may be made by dissolving the tunlametinib in the solvent of choice, which should be purged with an inert gas. Tunlametinib is soluble (≥10 mg/ml) in DMSO.

Description

Tunlametinib is a MEK1 inhibitor (IC₅₀ = 1.9 nM).¹ It is selective for MEK1 over a panel of other kinases at 10 μM. Tunlametinib selectively reduces proliferation in a panel of cancer cell lines expressing mutant K-Ras or B-RAF (IC₅₀s = 0.67-59.89 nM) over H1975 cells expressing wild-type K-Ras and B-RAF (IC₅₀ = >1,000 nM). It also inhibits the phosphorylation of ERK in A375 melanoma cells (IC₅₀ = 1.16 nM). *In vivo*, tunlametinib (1, 3, and 9 mg/kg) inhibits intratumoral ERK phosphorylation in an A375 mouse xenograft model and reduces tumor growth in the same model when administered at doses of 3 and 9 mg/kg.

Reference

1. Liu, Y., Cheng, Y., Huang, G., *et al.* Preclinical characterization of tunlametinib, a novel, potent, and selective MEK inhibitor. *Front. Pharmacol.* **14**, 1271268 (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, 02/16/2026

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