

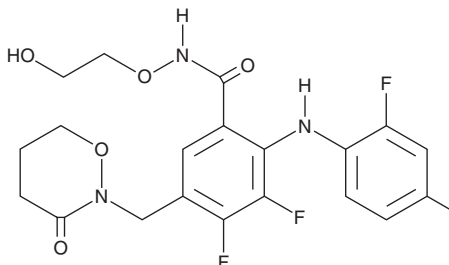
PRODUCT INFORMATION



RO4987655
Item No. 19488

CAS Registry No.: 874101-00-5
Formal Name: 3,4-difluoro-2-[(2-fluoro-4-iodophenyl)amino]-N-(2-hydroxyethoxy)-5-[(tetrahydro-3-oxo-2H-1,2-oxazin-2-yl)methyl]-benzamide

Synonym: CH4987655
MF: C₂₀H₁₉F₃IN₃O₅
FW: 565.3
Purity: ≥98%
UV/Vis.: λ_{max}: 280 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

RO4987655 is supplied as a crystalline solid. A stock solution may be made by dissolving the RO4987655 in the solvent of choice, which should be purged with an inert gas. RO4987655 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of RO4987655 in these solvents is approximately 10 and 15 mg/ml, respectively.

RO4987655 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, RO4987655 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. RO4987655 has a solubility of approximately 0.16 mg/ml in a 1:5 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

RO4987655 is a MEK inhibitor (IC₅₀ = 5 nM).¹ It is selective for MEK over 400 kinases at 10 μM. It inhibits proliferation of COLO 205, HT-29, QG56, MIA PaCa-2, and C32 cells with IC₅₀ values of 0.86, 1.7, 9.5, 3.3, and 8.4 nM, respectively. RO4987655 reduces tumor growth in a variety of mouse xenograft models and inhibits the phosphorylation of ERK in tumor tissue in an HT-29 mouse xenograft model when administered at a dose of 6.25 mg/kg per day. It acts synergistically with the mTOR inhibitor everolimus (Item No. 11597) to reduce tumor volume in an HCT116 mouse xenograft model.

Reference

1. Isshiki, Y., Kohchi, Y., Iikura, H., *et al.* Design and synthesis of novel allosteric MEK inhibitor CH4987655 as an orally available anticancer agent. *Bioorg. Med. Chem. Lett.* **21(6)**, 1795-1801 (2011).

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, 12/19/2022

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