

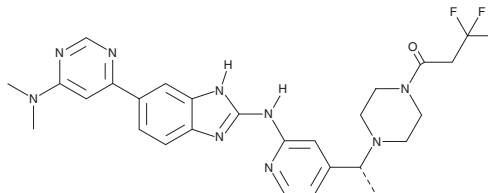
PRODUCT INFORMATION



BAY-985

Item No. 34234

CAS Registry No.: 2409479-29-2
Formal Name: 1-[4-[(1R)-1-[2-[[6-[6-(dimethylamino)-4-pyrimidinyl]-1H-benzimidazol-2-yl]amino]-4-pyridinyl]ethyl]-1-piperazinyl]-3,3,3-trifluoro-1-propanone
MF: C₂₇H₃₀F₃N₉O
FW: 553.6
Purity: ≥98%
UV/Vis.: λ_{max}: 266, 315 nm
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

BAY-985 is supplied as a solid. A stock solution may be made by dissolving the BAY-985 in the solvent of choice, which should be purged with an inert gas. BAY-985 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of BAY-985 in these solvents is approximately 25 and 30 mg/ml, respectively.

BAY-985 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, BAY-985 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. BAY-985 has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

BAY-985 is a dual inhibitor of TANK-binding kinase 1 (TBK1) and IκB kinase ε (IKKε; IC₅₀ = 2 nM for both).¹ It is selective for TBK1 and IKKε over FLT3, RSK4, DRAK1, and ULK1 (IC₅₀s = 123, 276, 311, and 7,390 nM, respectively). BAY-985 inhibits phosphorylation of interferon regulatory factor 3 (IRF3) in MDA-MB-231 cells expressing mouse IRF3 (IC₅₀ = 74 nM). It inhibits the proliferation of SK-MEL-2 cells *in vitro* (IC₅₀ = 900 nM) and reduces tumor weight in an SK-MEL-2 mouse xenograft model when administered at a dose of 200 mg/kg.

Reference

1. Lefranc, J., Schulze, V.K., Hillig, R.C., *et al.* Discovery of BAY-985, a highly selective TBK1/IKKε inhibitor. *J. Med. Chem.* **63**(2), 601-612 (2020).

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, 11/02/2022

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