

PRODUCT INFORMATION

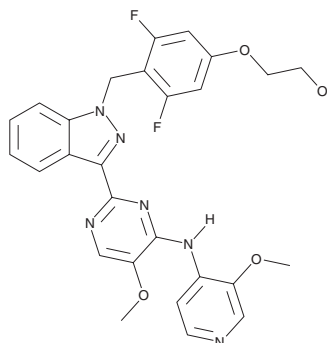


BAY-1816032

Item No. 35095

CAS Registry No.: 1891087-61-8
Formal Name: 2-[3,5-difluoro-4-[[3-[5-methoxy-4-[(3-methoxy-4-pyridinyl)amino]-2-pyrimidinyl]-1H-indazol-1-yl]methyl]phenoxy]-ethanol

MF: C₂₇H₂₄F₂N₆O₄
FW: 534.5
Purity: ≥98%
UV/Vis.: λ_{max}: 217, 257, 309 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-1816032 is supplied as a solid. A stock solution may be made by dissolving the BAY-1816032 in the solvent of choice, which should be purged with an inert gas. BAY-1816032 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of BAY-1816032 in these solvents is approximately 10 and 14 mg/ml, respectively. BAY-1816032 is slightly soluble in ethanol.

Description

BAY-1816032 is an inhibitor of the mitotic checkpoint serine/threonine protein kinase BUB1 (IC₅₀ = 6.1 nM).¹ It is selective for BUB1 over FLT3 (IC₅₀ = 4.2 μM) and 15 other kinases (IC₅₀s = >10 μM for all). BAY-1816032 inhibits proliferation of a variety of cancer cells (median IC₅₀ = 1.4 μM) and acts synergistically with the mitotic inhibitor paclitaxel (Item No. 10461) to increase the rate of chromosomal segregation defects during mitosis in HeLa cells. It reduces tumor growth in a SUM149 mouse xenograft model when administered at a dose of 25 mg/kg and has an additive effect on reducing tumor growth when used in combination with paclitaxel in the same model.

Reference

1. Siemeister, G., Mengel, A., Fernández-Montalván, A.E., *et al.* Inhibition of BUB1 kinase by BAY 1816032 sensitizes tumor cells toward taxanes, ATR, and PARP inhibitors *in vitro* and *in vivo*. *Clin. Cancer Res.* **25**(4), 1404-1414 (2019).

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, 12/02/2022

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