

PRODUCT INFORMATION



CB-6644

Item No. 36125

CAS Registry No.: 2316817-88-4
Formal Name: 5-chloro-N-[4-[[[6,7-dihydro-3-(methoxymethyl)-1-oxo-1H,5H-pyrazolo[1,2-a][1,2]benzodiazepin-2-yl]amino]-2,2-dimethyl-4-oxobutyl]-2-ethoxy-4-fluoro-benzamide

MF: C₂₉H₃₄ClFN₄O₅

FW: 573.1

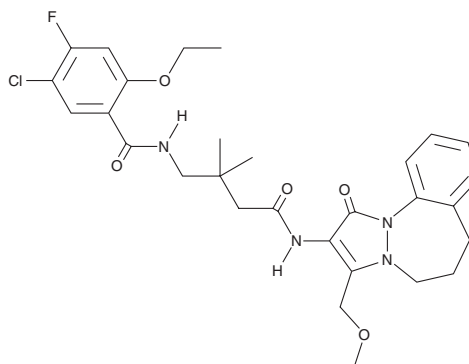
Purity: ≥98%

UV/Vis.: λ_{max}: 230 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

CB-6644 is supplied as a solid. A stock solution may be made by dissolving the CB-6644 in the solvent of choice, which should be purged with an inert gas. CB-6644 is soluble in DMSO.

Description

CB-6644 is an inhibitor of the RUVBL1/2 complex, a complex with ATPase activity that regulates several cellular processes, including chromatin remodeling and gene expression.¹ It inhibits the ATPase activity of the RUVBL1/2 complex (IC₅₀ = 15 nM) and is selective for the RUVBL1/2 complex over a panel of 224 ATPases at 10 μM. CB-6644 reduces viability in a panel of cancer cell lines with EC₅₀ values ranging from 41 to 785 nM. It induces apoptosis in HCT116 colon cancer cells when used at a concentration of 0.4 μM. CB-6644 (150 mg/kg) reduces tumor growth in Ramos Burkitt's lymphoma and RPMI-8226 multiple myeloma xenograft mouse models.

Reference

1. Assimon, V.A., Tang, Y., Vargas, J.D., *et al.* CB-6644 is a selective inhibitor of the RUVBL1/2 complex with anticancer activity. *ACS Chem Biol.* **14**(2), 236-244 (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, 10/25/2022

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