

PRODUCT INFORMATION



CCT369260

Item No. 41314

CAS Registry No.: 2253878-44-1

Formal Name: *rel*-5-[[5-chloro-2-[(3*R*,5*S*)-4,4-difluoro-3,5-dimethyl-1-piperidinyl]-4-pyrimidinyl]amino]-1,3-dihydro-3-(3-hydroxy-3-methylbutyl)-1-methyl-2*H*-benzimidazol-2-one

MF: C₂₄H₃₁ClF₂N₆O₂

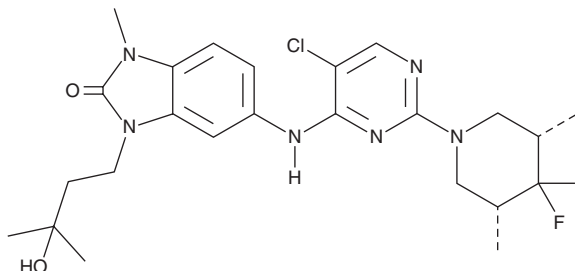
FW: 509.0

Purity: ≥95%

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

CCT369260 is supplied as a solid. A stock solution may be made by dissolving the CCT369260 in the solvent of choice, which should be purged with an inert gas. CCT369260 is sparingly soluble (1-10 mg/ml) in DMSO.

Description

CCT369260 is a degrader of Bcl-6.¹ It induces degradation of Bcl-6 in SU-DHL-4, OCI-LY1, and Karpas-422 cells with half-maximal degradation concentration (DC₅₀) values of 0.09, 0.049, and 0.062 μM, respectively. CCT369260 inhibits the growth of SU-DHL-4, OCI-LY1, and OCI-Ly3 cancer cells (GI₅₀s = 0.092, 0.035, and 1.6 μM, respectively). It reduces intratumoral Bcl-6 levels in an OCI-LY1 lymphoma mouse xenograft model for up to ten hours after administration of a 15 mg/kg dose.

Reference

1. Bellenie, B.R., Cheung, K.-M.J., Varela, A., *et al.* Achieving *in vivo* target depletion through the discovery and optimization of benzimidazolone BCL6 degraders. *J. Med. Chem.* **63**(8), 4047-4068 (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, 07/31/2024

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