

PRODUCT INFORMATION

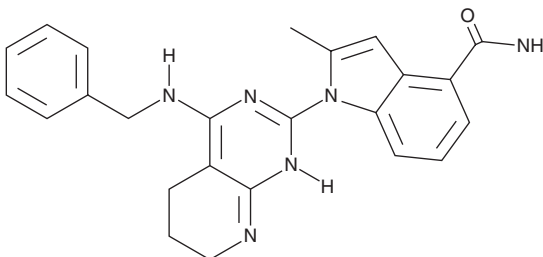


CB-5339

Item No. 46233

CAS Registry No.: 1863952-15-1
Formal Name: 2-methyl-1-[[5,6,7,8-tetrahydro-4-[(phenylmethyl)amino]pyrido[2,3-d]pyrimidin-2-yl]-1H-indole-4-carboxamide

MF: C₂₄H₂₄N₆O
FW: 412.5
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

CB-5339 is supplied as a solid. A stock solution may be made by dissolving the CB-5339 in the solvent of choice, which should be purged with an inert gas. CB-5339 is slightly soluble (0.1-1 mg/ml) in DMSO.

Description

CB-5339 is an inhibitor of the AAA-ATPase p97, also known as valosin-containing protein (VCP; IC₅₀ = 9 nM), and a derivative of CB-5083 (Item No. 19311).¹ It inhibits phosphodiesterase 6C (PDE6C) and accumulates in rat retina to a lesser extent than CB-5083. CB-5339 inhibits the growth of 16 primary acute myeloid leukemia (AML) cell lines (median IC₅₀ = 375 nM) and induces polyubiquitin protein accumulation and endoplasmic reticulum (ER) stress in MV4-11 AML cells. It decreases tumor cell bone marrow infiltration and increases survival in an MLL-AF9-driven patient-derived xenograft (PDX) mouse model of AML at 90 mg/kg.

Reference

1. Roux, B., Vaganay, C., Vargas, J.D., *et al.* Targeting acute myeloid leukemia dependency on VCP-mediated DNA repair through a selective second-generation small-molecule inhibitor. *Sci. Transl. Med.* **13**(587), eabg1168 (2021).

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, 08/31/2026

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