

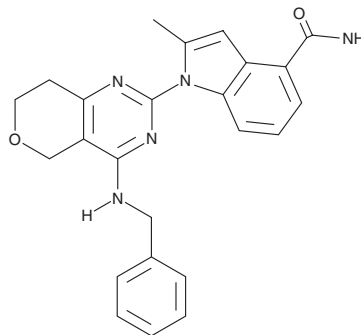
PRODUCT INFORMATION



CB-5083

Item No. 19311

CAS Registry No.: 1542705-92-9
Formal Name: 1-[7,8-dihydro-4-[(phenylmethyl)amino]-5H-pyrano[4,3-d]pyrimidin-2-yl]-2-methyl-1H-indole-4-carboxamide
MF: C₂₄H₂₃N₅O₂
FW: 413.5
Purity: ≥95%
UV/Vis.: λ_{max}: 245, 301 nm
Supplied as: A crystalline 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-5083 is supplied as a crystalline solid. A stock solution may be made by dissolving the CB-5083 in the solvent of choice, which should be purged with an inert gas. CB-5083 is sparingly soluble in warmed ethanol.

Description

CB-5083 is an orally bioavailable inhibitor of p97 (IC₅₀ = 11 nM), a transitional endoplasmic reticulum ATPase also known as valosin-containing protein (VCP).^{1,2} It disrupts cellular protein homeostasis and demonstrates anti-tumor activity in hematological and solid tumor models. CB-5083 causes accumulation of poly-ubiquitinated proteins, retention of endoplasmic reticulum-associated degradation (ERAD) substrates, and activation of the apoptotic arm of the unfolded protein response.^{1,2}

References

1. Anderson, D.J., Le Moigne, R., Djjakovic, S., *et al.* Targeting the AAA ATPase p97 as an approach to treat cancer through disruption of protein homeostasis. *Cancer Cell*. **28(5)**, 653-665 (2015).
2. Zhou, H.-J., Wang, J., Yao, B., *et al.* Discovery of a first-in-class, potent, selective, and orally bioavailable inhibitor of the p97 AAA ATPase (CB-5083). *J. Med. Chem.* **58(24)**, 9480-9497 (2015).

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, 11/22/2022

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