

PRODUCT INFORMATION



RP 1664

Item No. 44283

CAS Registry No.: 2980682-00-4
Formal Name: 5-cyclopropyl-N²-(2,6-difluoro-4-(methylsulfonyl)phenyl)-N²-methyl-6-(1-methyl-1H-imidazol-4-yl)-N⁴-(5-methyl-1H-pyrazol-3-yl)pyrimidine-2,4-diamine

MF: C₂₃H₂₄F₂N₈O₂S

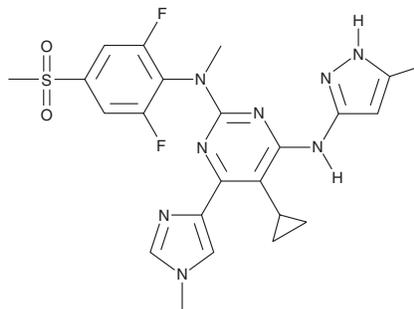
FW: 514.6

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

RP 1664 is supplied as a solid. A stock solution may be made by dissolving the RP 1664 in the solvent of choice, which should be purged with an inert gas. RP 1664 is sparingly slightly soluble (0.1-1 mg/ml) in acetonitrile and DMSO.

Description

RP 1664 is an inhibitor of polo-like kinase 4 (Plk4; IC₅₀ = 3 nM).¹ It induces centriole depletion in hTERT RPE-1 retinal pigment cells (EC₅₀ = 0.13 μM). RP 1664 induces cytotoxicity in MCF-7 breast cancer cells (EC₅₀ = 0.051 μM). It reduces tumor volume in a CAL-148 breast cancer mouse xenograft model when administered at doses of 11 or 21 mg/kg twice per day.

Reference

- Vallée, F., Casás-Selves, M., Bubenik, M., *et al.* Discovery of RP-1664: A first-in-class orally bioavailable, selective PLK4 inhibitor. *J. Med. Chem.* **68(11)**, 10631-10647 (2025).

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, 09/22/2025

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