

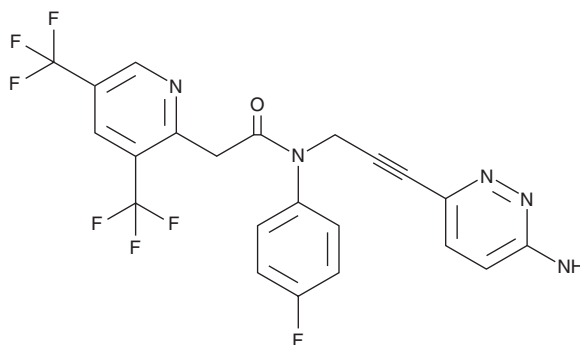
PRODUCT INFORMATION



RP 6685

Item No. 42127

CAS Registry No.: 2832047-80-8
Formal Name: N-[3-(6-amino-3-pyridazinyl)-2-propyn-1-yl]-N-(4-fluorophenyl)-3,5-bis(trifluoromethyl)-2-pyridineacetamide
MF: C₂₂H₁₄F₇N₅O
FW: 497.4
Purity: ≥98%
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 6685 is supplied as a solid. A stock solution may be made by dissolving the RP 6685 in the solvent of choice, which should be purged with an inert gas. RP 6685 is soluble (≥10 mg/ml) in DMSO.

Description

RP 6685 is an inhibitor of DNA polymerase θ (IC_{50} = 0.55 nM).¹ It is selective for DNA polymerase θ over DNA polymerase α , $-\epsilon$, $-\gamma$, and $-\nu$ (IC_{50} s = >100 μ M for all). RP 6685 inhibits double-strand break repair in adeno-associated virus 1 (AAVS1) loci in HCT116 cells (IC_{50} = 0.5 μ M) and the proliferation of HCT116 BRCA2^{-/-} cells (IC_{50} = 0.32 μ M). *In vivo*, RP 6685 (80 mg/kg) reduces tumor volume in an HCT116 BRCA2^{-/-} mouse xenograft model.

Reference

1. Bubenik, M., Mader, P., Mochirian, P., *et al.* Identification of RP-6685, an orally bioavailable compound that inhibits the DNA polymerase activity of Pol θ . *J. Med. Chem.* **65**(19), 13198-13215 (2022).

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, 10/30/2024

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