

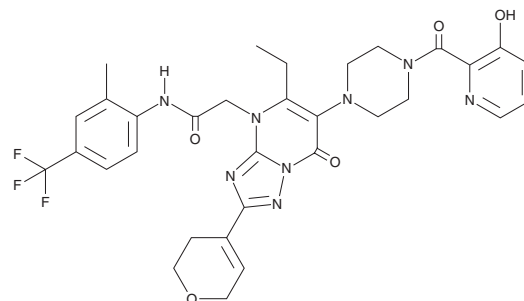
PRODUCT INFORMATION



Werner Syndrome RecQ Helicase-IN-4

Item No. 41679

CAS Registry No.: 2869954-53-8
Formal Name: 2-(3,6-dihydro-2H-pyran-4-yl)-5-ethyl-6-[4-[(3-hydroxy-2-pyridinyl)carbonyl]-1-piperazinyl]-N-[2-methyl-4-(trifluoromethyl)phenyl]-7-oxo-[1,2,4]triazolo[1,5-a]pyrimidine-4(7H)-acetamide
Synonyms: Werner Syndrome RecQ Helicase Inhibitor 4, WS RecQ Helicase-IN-4, WS RecQ Helicase Inhibitor 4
MF: C₃₂H₃₃F₃N₈O₅
FW: 666.7
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

Werner syndrome RecQ helicase-IN-4 is supplied as a solid. A stock solution may be made by dissolving the werner syndrome RecQ helicase-IN-4 in the solvent of choice, which should be purged with an inert gas. Werner syndrome RecQ helicase-IN-4 is sparingly soluble (1-10 mg/ml) in DMSO.

Description

Werner syndrome RecQ helicase-IN-2 is an inhibitor of the DNA helicase Werner syndrome protein (WRN; IC₅₀ = 60 nM).¹ It selectively inhibits the growth of SW48 colorectal cancer cells over DLD-1 colorectal cancer cells lacking WRN (DLD-1-WRN-KO; GI₅₀s = 70 and >10,000 nM, respectively). Werner syndrome RecQ helicase-IN-2 (240 mg/kg per day) reduces tumor volume in an SW48 mouse xenograft model.

Reference

1. Bordas, V., Brun, J., Decker, A., *et al.* Triazolo-pyrimidine analogues for treating diseases connected to the inhibitor of Werner syndrome RecQ helicase (WRN). *Novartis AG, WO2022/249060AI* (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, 10/31/2024

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