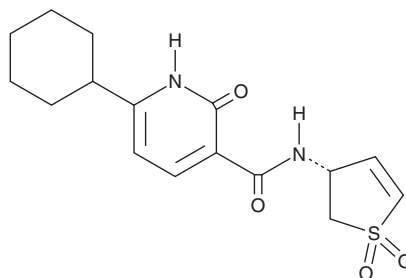


PRODUCT INFORMATION



GSK_WRN4
Item No. 45316

CAS Registry No.: 2923008-66-4
Formal Name: 6-cyclohexyl-N-[(3R)-2,3-dihydro-1,1-dioxido-3-thienyl]-1,2-dihydro-2-oxo-3-pyridinecarboxamide
MF: C₁₆H₂₀N₂O₄S
FW: 336.4
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

GSK_WRN4 is supplied as a solid. A stock solution may be made by dissolving the GSK_WRN4 in the solvent of choice, which should be purged with an inert gas. GSK_WRN4 is slightly soluble (0.1-1 mg/ml) in acetonitrile and DMSO.

Description

GSK_WRN4 is a covalent inhibitor of the DNA helicase Werner syndrome protein (WRN; IC₅₀ = 25 nM).¹ It inhibits tumor growth and increases intratumoral levels of the DNA damage markers p21, phosphorylated γH2AX, and phosphorylated KAP-1 in an SW48 colorectal cancer mouse xenograft model when administered at a dose of 300 mg/kg.

Reference

1. Picco, G., Rao, Y., Al Saedi, A., *et al.* Novel WRN helicase inhibitors selectively target microsatellite-unstable cancer cells. *Cancer Discov.* **14**(8), 1457-1475 (2024).

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, 03/18/2026

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