

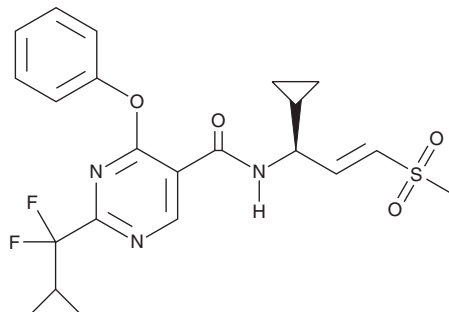
PRODUCT INFORMATION



WRN Inhibitor 8

Item No. 43724

CAS Registry No.: 3026502-95-1
Formal Name: 2-(cyclopropyl)difluoromethyl)-N-[(1S,2E)-1-cyclopropyl-3-(methylsulfonyl)-2-propen-1-yl]-4-phenoxy-5-pyrimidinecarboxamide
MF: C₂₂H₂₃F₂N₃O₄S
FW: 463.5
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

WRN inhibitor 8 is supplied as a solid. A stock solution may be made by dissolving the WRN inhibitor 8 in the solvent of choice, which should be purged with an inert gas. WRN inhibitor 8 is sparingly soluble (1-10 mg/ml) in DMSO.

Description

WRN inhibitor 8 is an inhibitor of the DNA helicase Werner syndrome protein (WRN; IC₅₀ = 48.3 nM).¹ It inhibits the growth of HCT116 and SW480 colorectal cancer cells (GI₅₀s = 0.01699 and 13.64 μM, respectively).

Reference

1. Kikuchi, S., Lam, B., Green, J., *et al.* Pharmaceutical compositions comprising WRN helicase inhibitors. *Vividion Therapeutics Inc. WO2024010782A1* (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, 07/29/2025

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