

PRODUCT INFORMATION



WDR5-MYC Probe 12

Item No. 39304

Formal Name: 3',5'-difluoro-4'-((6-methyl-4-oxo-1-phenyl-1,4-dihydropyridazine-3-carboxamido)methyl)-[1,1'-biphenyl]-3-sulfonyl fluoride

MF: C₂₅H₁₈F₃N₃O₄S

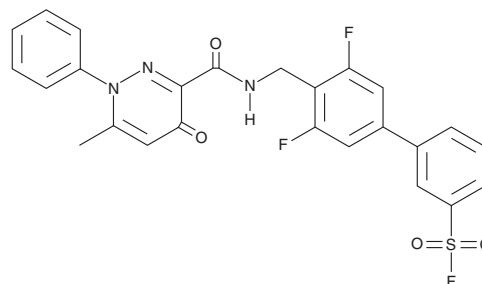
FW: 513.5

Purity: ≥98%

Supplied as: A solid

Storage: -20°C

Stability: ≥3 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

WDR5-MYC probe 12 is supplied as a solid. A stock solution may be made by dissolving the WDR5-MYC probe 12 in the solvent of choice, which should be purged with an inert gas. WDR5-MYC probe 12 is soluble in acetonitrile and DMSO.

Description

WDR5-MYC probe 12 is an inhibitor of the protein-protein interaction between the master transcription factor Myc and WD repeat domain 5 (WDR5; IC₅₀ = 0.39 μM).¹ It reduces mRNA expression of the Myc and WDR5 co-binding target genes ribosomal protein L17 (*RPL17*), *RPL23*, *RPL35*, and *RPL37A* in HL-60 cells. WDR5-MYC probe 12 also inhibits the growth of HL-60, NB4, Ramos (RA 1), COR-L279, and NCI H520 cancer cells (GI₅₀s = 4.2-9.4 μM).

Reference

1. Ding, J., Liu, L., Chiang, Y.L., *et al.* Discovery and structure-based design of inhibitors of the WD repeat-containing protein 5 (WDR5)-MYC interaction. *J. Med. Chem.* **66**(12), 8310-8323 (2023).

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, 12/22/2023

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