

PRODUCT INFORMATION

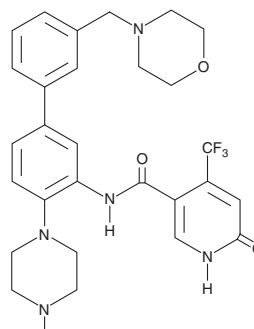


OICR-9429

Item No. 16095

CAS Registry No.: 1801787-56-3
Formal Name: N-(4-(4-methylpiperazin-1-yl)-3'-(morpholinomethyl)-[1,1'-biphenyl]-3-yl)-6-oxo-4-(trifluoromethyl)-1,6-dihydropyridine-3-carboxamide

MF: C₂₉H₃₂F₃N₅O₃
FW: 555.6
Purity: ≥98%
UV/Vis.: λ_{max}: 259 nm
Supplied as: A crystalline 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

OICR-9429 is supplied as a crystalline solid. A stock solution may be made by dissolving the OICR-9429 in the solvent of choice, which should be purged with an inert gas. OICR-9429 is soluble in organic solvents such as DMSO and methanol. The solubility of OICR-9429 in these solvents is approximately 0.5 mg/ml (with gentle warming).

OICR-9429 is sparingly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

WD-repeat protein 5 (WDR5) is a scaffold protein commonly involved in the formation of nucleosome-modifying protein complexes with histones.¹ For example, WDR5 is a component of a mixed-lineage leukemia (MLL) methyltransferase complex that targets histone 3.² WDR5 can also be a component of histone acetyltransferase complexes and can directly bind methylated as well as unmodified histones.^{1,3} OICR-9429 is a chemical probe that inhibits the interaction of WDR5 with peptide regions of MLL and Histone 3, demonstrating >100-fold selectivity over other chromatin reader domains, methyltransferases, and other non-epigenetic targets. It selectively binds to WDR5 (K_d = 24-52 nM) and disrupts its interaction with MLL1 and RbBP5 in cells (IC₅₀ = < 1 μM). For more information on OICR-9429 please visit the Structural Genomics Consortium (SGC). The negative control, OICR-0547, for OICR-9429 is also available exclusively through the SGC.

References

1. Migliori, V., Mapelli, M., and Guccione, E. On WD40 proteins: Propelling our knowledge of transcriptional control? *Epigenetics* **7**(8), 815-822 (2012).
2. Guccione, E., Bassi, C., Casadio, F., et al. Methylation of histone H3R2 by PRMT6 and H3K4 by an MLL complex are mutually exclusive. *Nature* **449**, 933-937 (2007).
3. Guelman, S., Kozuka, K., Mao, Y., et al. The double-histone-acetyltransferase complex ATAC is essential for mammalian development. *Mol. Cell. Biol.* **29**(5), 1176-1188 (2009).

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, 01/03/2023

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