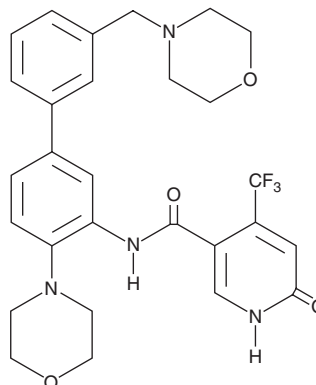


PRODUCT INFORMATION

OICR-0547

Item No. 18180

CAS Registry No.: 1801873-49-3
Formal Name: 1,6-dihydro-N-[4-(4-morpholinyl)-3'-(4-morpholinylmethyl)[1,1'-biphenyl]-3-yl]-6-oxo-4-(trifluoromethyl)-3-pyridinecarboxamide
MF: C₂₈H₂₉F₃N₄O₄
FW: 542.6
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

OICR-0547 is supplied as a solid. A stock solution may be made by dissolving the OICR-0547 in the solvent of choice, which should be purged with an inert gas. OICR-0547 is sparingly soluble (1-10 mg/ml) in DMSO.

Description

OICR-0547 is an inactive derivative of OICR-9429 (Item No. 16095), an inhibitor of the protein-protein interaction between WD40 repeat protein 5 (WDR5) and mixed-lineage leukemia (MLL).¹ It has been used as a negative control for the protein-protein inhibition activity of OICR-9429. See the Structural Genomics Consortium (SGC) website for more information.

Reference

1. Grebien, F., Vedadi, M., Getlik, M., *et al.* Pharmacological targeting of the Wdr5-MLL interaction in C/EBPα N-terminal leukemia. *Nat. Chem. Biol.* (2015).

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, 02/20/2025

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