

PRODUCT INFORMATION

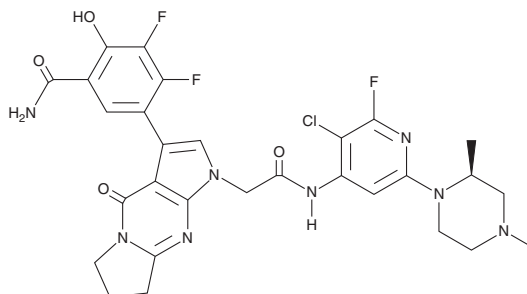


OICR-12694

Item No. 44288

CAS Registry No.: 2360625-97-2
Formal Name: 3-[5-(aminocarbonyl)-2,3-difluoro-4-hydroxyphenyl]-N-[3-chloro-6-[(2S)-2,4-dimethyl-1-piperazinyl]-2-fluoro-4-pyridinyl]-4,6,7,8-tetrahydro-4-oxo-1H-dipyrrolo[1,2-a:2',3'-d]pyrimidine-1-acetamide

Synonym: JNJ-65234637
MF: C₂₉H₂₈ClF₃N₈O₄
FW: 645.0
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

OICR-12694 is supplied as a solid. A stock solution may be made by dissolving the OICR-12694 in the solvent of choice, which should be purged with an inert gas. OICR-12694 is sparingly soluble (1-10 mg/ml) in DMSO.

Description

OICR-12694 is an inhibitor of Bcl-6.¹ It binds to the BTB domain on Bcl-6 ($K_d = 5$ nM in a surface plasmon resonance (SPR) assay). OICR-12694 decreases transcription of Bcl-6-targeted genes in a reporter assay using SU-DHL-4 cells ($EC_{50} = 89$ nM). It reduces the proliferation of Karpas-422 large B cell lymphoma cells ($IC_{50} = 92$ nM).

Reference

1. Mamai, A., Chau, A.M., Wilson, B.J., *et al.* Discovery of OICR12694: A novel, potent, selective, and orally bioavailable BCL6 BTB inhibitor. *ACS Med. Chem. Lett.* **14**(2), 199-210 (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, 08/19/2025

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