

PRODUCT INFORMATION

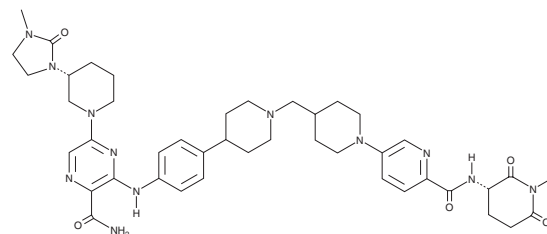


NX-5948

Item No. 43484

CAS Registry No.: 2649400-34-8
Formal Name: 3-[[4-[1-[[1-[6-[[[(3S)-2,6-dioxo-3-piperidinyl]amino]carbonyl]-3-pyridinyl]-4-piperidinyl]methyl]-4-piperidinyl]phenyl]amino]-5-[(3R)-3-(3-methyl-2-oxo-1-imidazolidiny)]-1-piperidinyl]-2-pyrazinecarboxamide

MF: C₄₂H₅₄N₁₂O₅
FW: 807.0
Purity: ≥98%
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

NX-5948 is supplied as a crystalline solid. A stock solution may be made by dissolving the NX-5948 in the solvent of choice, which should be purged with an inert gas. NX-5948 is sparingly soluble (1-10 mg/ml) in DMSO.

Description

NX-5948 is a proteolysis-targeting chimera (PROTAC) that drives the degradation of Bruton's tyrosine kinase (BTK).^{1,2} It induces the degradation of BTK in OCI-LY10 diffuse large B cell lymphoma (DLBCL) cells with a half-maximal degradation concentration (DC₅₀) value of 0.07 nM.² NX-5948 inhibits the proliferation of wild-type BTK-expressing OCI-LY10 and TMD8 DLBCL cells and mutant BTK^{C481S}-expressing TMD8 cells (IC₅₀s = 0.1, 0.6, and 16 nM, respectively).

References

1. Wang, F., Sun, W., Ma, Z., *et al.* A validated liquid chromatography-tandem mass spectrometry assay for the determination of NX-5948 in beagle dog plasma: Application to a pharmacokinetic study. *Biomed. Chromatogr.* **39(3)**, e6090 (2025).
2. Ren, J., Huang, Y., Wang, J., *et al.* Discovery of TQ-3959 as a potent and orally bioavailable BTK PROTAC degrader incorporating a novel benzisoxazole-based CRBN ligand for the treatment of B-cell malignancies. *J. Med. Chem.* **Online ahead of print**, (2025).

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, 08/07/2025

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