

PRODUCT INFORMATION

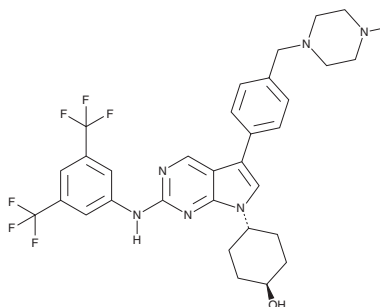


UNC9426

Item No. 44849

CAS Registry No.: 3051784-94-9
Formal Name: *trans*-4-[2-[[3,5-bis(trifluoromethyl)phenyl]amino]-5-[4-[(4-methyl-1-piperazinyl)methyl]phenyl]-7H-pyrrolo[2,3-d]pyrimidin-7-yl]-cyclohexanol

MF: C₃₂H₃₄F₆N₆O
FW: 632.7
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

UNC9426 is supplied as a solid. A stock solution may be made by dissolving the UNC9426 in the solvent of choice, which should be purged with an inert gas. UNC9426 is slightly soluble (0.1-1 mg/ml) in ethanol. UNC9426 is soluble (≥10 mg/ml) in DMSO.

Description

UNC9426 is an inhibitor of the TAM family kinase Tyro3 (IC₅₀ = 2.1 nM).¹ It is selective for Tyro3 over MerTK and Axl (IC₅₀s = 580 and 190 nM, respectively) and 74 other kinases in a panel at 1 μM. UNC9426 (10 μM) reduces ADP- or collagen-induced platelet aggregation in isolated human platelet-rich plasma. It does not increase total bleeding time in mice after tail tip transection when administered at a dose of 10 mg/kg. UNC9426 also reduces the uptake of phosphatidylserine-coated microparticles, a model of apoptotic cells, by THP-1 macrophages in an efferocytosis assay in a concentration-dependent manner and reduces colony formation in NCI H1650 or H1975 non-small cell lung cancer (NSCLC) cells (EC₅₀s = 170 and 197 nM, respectively).

Reference

1. Kong, D., Yang, X., Judd, S., *et al.* UNC9426, a potent and orally bioavailable TYRO3-specific inhibitor. *J. Med. Chem.* **68**(6), 665-6682 (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, 01/26/2026

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