

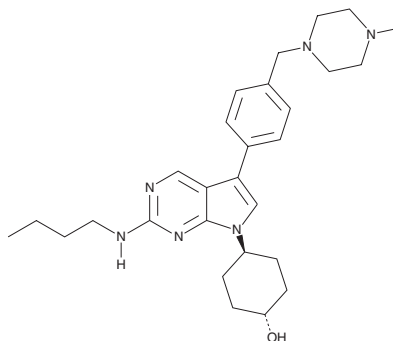
# PRODUCT INFORMATION



## UNC2025

Item No. 16613

**CAS Registry No.:** 1429881-91-3  
**Formal Name:** *trans*-4-[2-(butylamino)-5-[4-[(4-methyl-1-piperazinyl)methyl]phenyl]-7H-pyrrolo[2,3-d]pyrimidin-7-yl]-cyclohexanol  
**MF:** C<sub>28</sub>H<sub>40</sub>N<sub>6</sub>O  
**FW:** 476.7  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 222, 254, 290, 337 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

UNC2025 is supplied as a crystalline solid. A stock solution may be made by dissolving the UNC2025 in the solvent of choice, which should be purged with an inert gas. UNC2025 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of UNC2025 in ethanol and DMF is approximately 30 mg/ml, and approximately 10 mg/ml in DMSO.

UNC2025 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, UNC2025 should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. UNC2025 has a solubility of approximately 0.1 mg/ml in a 1:9 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

UNC2025 is a potent, orally bioavailable inhibitor of the tyrosine kinases Mer and Flt3 (IC<sub>50</sub>s = 0.74 and 0.80 nM, respectively).<sup>1</sup> It less potently inhibits Axl and Tyro3 (IC<sub>50</sub>s = 14 and 17 nM, respectively) and a panel of related kinases.<sup>1</sup> UNC2025 blocks colony formation of Mer- and Flt3-dependent tumor cell lines and inhibits Mer phosphorylation in bone marrow leukemia cells *in vivo*.<sup>1</sup> Pharmacokinetic studies suggest that orally administered UNC2025 at 3 mg/kg will suffice to provide 90% inhibition of Mer and Flt3 at 30 min.<sup>1</sup>

### References

1. Zhang, W., DeRyckere, D., Hunter, D., *et al.* UNC2025, a potent and orally bioavailable MER/FLT3 dual inhibitor. *J. Med. Chem.* **57**(16), 7031-7041 (2014).

#### 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, 12/08/2022

#### 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