

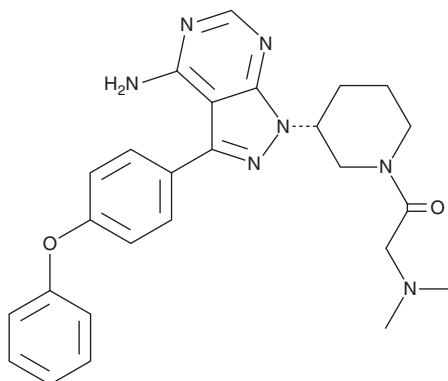
PRODUCT INFORMATION



CHMFL-FLT3-122

Item No. 43137

CAS Registry No.: 1839150-56-9
Formal Name: 1-[(3R)-3-[4-amino-3-(4-phenoxyphenyl)-1H-pyrazolo[3,4-d]pyrimidin-1-yl]-1-piperidinyl]-2-(dimethylamino)ethanone
Synonym: HYML-122
MF: C₂₆H₂₉N₇O₂
FW: 471.6
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

CHMFL-FLT3-122 is supplied as a solid. A stock solution may be made by dissolving the CHMFL-FLT3-122 in the solvent of choice, which should be purged with an inert gas. CHMFL-FLT3-122 is slightly soluble (0.1-1 mg/ml) in acetonitrile and DMSO.

Description

CHMFL-FLT3-122 is an inhibitor of FMS-related tyrosine kinase 3 (FLT3; IC₅₀ = 40 nM).¹ It is selective for FLT3 over Bruton's tyrosine kinase (BTK) and C-Kit (IC₅₀s = 421 and 559 nM, respectively) and a panel of 468 additional kinases at 1 μM. CHMFL-FLT3-122 inhibits the growth of MV4-11, MOLM-13, and MOLM-14 acute myeloid leukemia (AML) cells containing FLT3 bearing an internal-tandem duplication (FLT3-ITD; GI₅₀s = 22, 21, and 42 nM, respectively). *In vivo*, CHMFL-FLT3-122 (50 mg/kg) inhibits tumor growth in an MV4-11 mouse xenograft model.

Reference

1. Li, X., Wang, A., Yu, K., *et al.* Discovery of (R)-1-(3-(4-amino-3-(4-phenoxyphenyl)-1H-pyrazolo[3,4-d]pyrimidin-1-yl)piperidin-1-yl)-2-(dimethylamino)ethanone (CHMFL-FLT3-122) as a potent and orally available FLT3 kinase inhibitor for FLT3-ITD positive acute myeloid leukemia. *J. Med. Chem.* **58**(24), 9625-9638 (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, 04/08/2025

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