

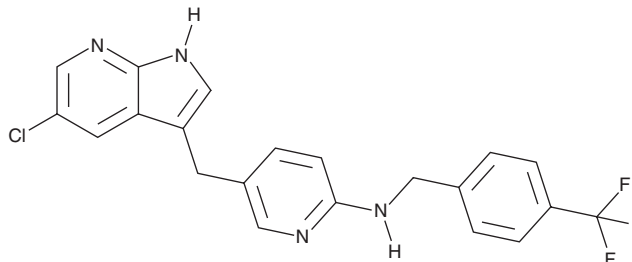
PRODUCT INFORMATION



FLT3-IN-2

Item No. 30070

CAS Registry No.: 923562-23-6
Formal Name: 5-[(5-chloro-1H-pyrrolo[2,3-b]pyridin-3-yl)methyl]-N-[[4-(trifluoromethyl)phenyl]methyl]-2-pyridinamine
MF: C₂₁H₁₆ClF₃N₄
FW: 416.8
Purity: ≥98%
UV/Vis.: λ_{max}: 235, 305 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

FLT3-IN-2 is supplied as a solid. Aqueous solutions of FLT3-IN-2 can be prepared by directly dissolving the solid in aqueous buffers. The solubility of FLT3-IN-2 in PBS, pH 7.2, is approximately 0.25 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

FLT3-IN-2 is an inhibitor of FMS-related tyrosine kinase 3 (FLT3), c-Kit, and c-FMS (IC₅₀s = <1 μM for all).^{1,2}

References

1. Zhang, C., Zhang, J., Ibrahim, P., *et al.* Compounds modulating c-fms and/or c-kit activity and uses therefor. *Plexxikon, Inc. WO2008/064265A2* (2008).
2. Zhang, C., Bollag, G., Habets, G., *et al.* Kinase modulation and indications therefor. *Plexxikon, Inc. US2013/0274259A1* (2013).

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, 05/20/2020

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