

PRODUCT INFORMATION



B-RAF IN 1

Item No. 23434

CAS Registry No.: 950736-05-7
Formal Name: N-[3-[3-[4-[(dimethylamino)methyl]phenyl]pyrazolo[1,5-a]pyrimidin-7-yl]phenyl]-3-(trifluoromethyl)-benzamide

Synonym: B-RAF Inhibitor 1

MF: C₂₉H₂₄F₃N₅O

FW: 515.5

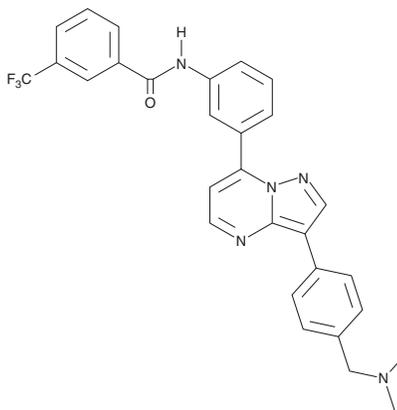
Purity: ≥98%

UV/Vis.: λ_{max}: 269 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

B-RAF IN 1 is supplied as a crystalline solid. A stock solution may be made by dissolving the B-RAF IN 1 in the solvent of choice. B-RAF IN 1 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of B-RAF IN 1 in these solvents is approximately 1, 33, and 14 mg/ml, respectively.

Description

B-RAF IN 1 is an inhibitor of B-RAF (IC₅₀ = 24 nM) that also inhibits C-Raf (IC₅₀ = 25 nM).¹ It is selective over 13 other kinases, including PKCα, IKKβ, and PI3Kα, at concentrations greater than 2 μM, but does not inhibit p38α and CAMKII (IC₅₀s = 216 and 822 nM, respectively). B-RAF IN 1 binds to B-RAF in the inactive conformation based on co-crystallization with the wild-type enzyme. It inhibits proliferation of WM 266-4 and HT29 cells with IC₅₀ values of 920 and 780 nM, respectively.

Reference

- Berger, D.M., Torres, N., Dutia, M., *et al.* Non-hinge-binding pyrazolo[1,5-a]pyrimidines as potent B-Raf kinase inhibitors. *Bioorg. Med. Chem. Lett.* **19(23)**, 6519-6523 (2009).

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, 12/06/2022

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