

PRODUCT INFORMATION

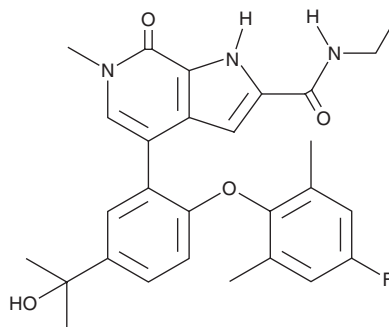


ABBV-744

Item No. 30470

CAS Registry No.: 2138861-99-9
Formal Name: N-ethyl-4-[2-(4-fluoro-2,6-dimethylphenoxy)-5-(1-hydroxy-1-methylethyl)phenyl]-6-methyl-7-oxo-1H-pyrrolo[2,3-c]pyridine-2-carboxamide

MF: C₂₈H₃₀FN₃O₄
FW: 491.6
Purity: ≥98%
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

ABBV-744 is supplied as a crystalline solid. A stock solution may be made by dissolving the ABBV-744 in the solvent of choice, which should be purged with an inert gas. ABBV-744 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of ABBV-744 in these solvents is approximately 30 mg/ml.

ABBV-744 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, ABBV-744 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. ABBV-744 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

ABBV-744 is an inhibitor of the bromodomain and extra-terminal domain (BET) family protein bromodomain 2 (BD2; K_ds = 4.6, 4.9, 1.6, and 1 nM for BRD2, -3, -4, and -T, respectively).¹ It is selective for BD2 over BD1 bromodomains in a time-resolved FRET (TR-FRET) assay (K_ds = 1,162, 3,140, 521, and 917 nM for BRD2, -3, -4, and -T, respectively). ABBV-744 inhibits proliferation of SKM-1 leukemia cells (IC₅₀ = 6.7 nM). It induces cell cycle arrest at the G₁ phase in LNCaP prostate cancer cells.² ABBV-744 (18.8 mg/kg) reduces tumor volume in a SKM-1 mouse xenograft model.¹ It also reduces tumor growth in LNCaP and MDA PCa 2b mouse xenograft models when administered at doses of 1 and 4.7 mg/kg.²

References

1. Sheppard, G.S., Wang, L., Fidanze, S.D., *et al.* Discovery of N-Ethyl-4-[2-(4-fluoro-2,6-dimethyl-phenoxy)-5-(1-hydroxy-1-methyl-ethyl)phenyl]-6-methyl-7-oxo-1H-pyrrolo[2,3-c]pyridine-2-carboxamide (ABBV-744), a BET bromodomain inhibitor with selectivity for the second bromodomain. *J. Med. Chem.* **63**(10), 5585-5623 (2020).
2. Faivre, E., McDaniel, K.F., Albert, D.H., *et al.* Selective inhibition of the BD2 bromodomain of BET proteins in prostate cancer. *Nature.* **578**(7994), 306-310 (2020).

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, 12/12/2022

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