

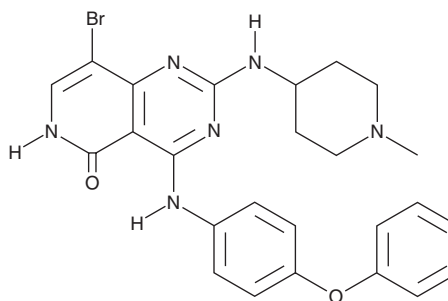
PRODUCT INFORMATION



G-749

Item No. 22139

CAS Registry No.: 1457983-28-6
Formal Name: 8-bromo-2-[(1-methyl-4-piperidinyl)amino]-4-[(4-phenoxyphenyl)amino]-pyrido[4,3-d]pyrimidin-5(6H)-one
MF: C₂₅H₂₅BrN₆O₂
FW: 521.4
Purity: ≥98%
UV/Vis.: λ_{max}: 261 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

G-749 is supplied as a crystalline solid. A stock solution may be made by dissolving the G-749 in the solvent of choice, which should be purged with an inert gas. G-749 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of G-749 in these solvents is approximately 3 and 16 mg/ml, respectively.

G-749 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, G-749 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. G-749 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

G-749 is an inhibitor of Fms-like tyrosine receptor kinase 3 (FLT3) with IC₅₀ values ranging from 2.1 to 38.1 nM for wild-type and mutant (constitutively active) FLT3s in Ba/F3 cells expressing recombinant receptors.¹ It also inhibits growth of human leukemia cell lines that express FLT3-ITD mutant kinase (IC₅₀s = 3.5 and 7.5 nM for MV4-11 and Molm-14 cells, respectively) but has no effect on cells that lack FLT3 expression or that express wild-type FLT3. G-749 reduces tumor growth and increases survival in a dose-dependent manner in MV4-11 and Molm-14 murine leukemia xenograft models. It reduces growth of bone marrow blasts derived from acute myeloid leukemia patients, showing more potent inhibition of blasts expressing the FLT3-ITD mutant receptor.

Reference

1. Lee, H.K., Kim, H.W., Lee, I.Y., et al. G-749, a novel FLT3 kinase inhibitor, can overcome drug resistance for the treatment of acute myeloid leukemia. *Blood* **123**(14), 2209-2219 (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, 11/14/2022

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