

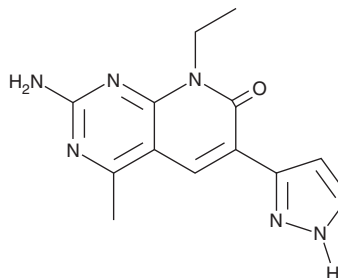
PRODUCT INFORMATION



Voxtalib

Item No. 30323

CAS Registry No.: 934493-76-2
Formal Name: 2-amino-8-ethyl-4-methyl-6-(1H-pyrazol-3-yl)-pyrido[2,3-d]pyrimidin-7(8H)-one
Synonyms: SAR245409, XL765
MF: C₁₃H₁₄N₆O
FW: 270.3
Purity: ≥98%
UV/Vis.: λ_{max}: 220, 362 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

Voxtalib is supplied as a crystalline solid. A stock solution may be made by dissolving the voxtalib in the solvent of choice, which should be purged with an inert gas. Voxtalib is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of voxtalib in these solvents is approximately 10 and 1 mg/ml, respectively.

Voxtalib is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, voxtalib should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Voxtalib has a solubility of approximately 0.16 mg/ml in a 1:5 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Voxtalib is a dual inhibitor of PI3K and mammalian target of rapamycin complex (mTORC; IC₅₀s = 39, 110, 43, 9, 160, and 910 nM for PI3Kα, PI3Kβ, PI3Kδ, PI3Kγ, mTORC1, and mTORC2, respectively).¹ It is selective for these kinases over vacuolar protein sorting 34 (VPS34; IC₅₀ = 9,100 nM) and a panel of 130 additional protein kinases at 1.5 μM but does inhibit DNA protein kinase (DNA-PK; IC₅₀ = 150 nM). Voxtalib decreases EGF-induced production of PIP₃, a product of PI3K-mediated PIP₂ metabolism, in PC3 prostate cancer cells with an IC₅₀ value of 290 nM. It also reduces phosphorylation of the PI3K and mTORC targets Akt and ribosomal protein S6 (S6RP) in the same model (IC₅₀s = 250 and 120 nM, respectively). Voxtalib inhibits proliferation and colony formation of PC3 cells (IC₅₀s = 1,800 and 270 nM, respectively). It reduces tumor growth and increases survival in a GBM-39 glioblastoma multiforme (GBM) mouse xenograft model when administered at a dose of 30 mg/kg twice per day alone or in combination with temozolomide (Item No. 14163).²

References

1. Yu, P., Laird, A.D., Du, X., *et al.* Characterization of the activity of the PI3K/mTOR inhibitor XL765 (SAR245409) in tumor models with diverse genetic alterations affecting the PI3K pathway. *Mol. Cancer Ther.* **13**(5), 1078-1091 (2014).
2. Prasad, G., Sottero, T., Yang, X., *et al.* Inhibition of PI3K/mTOR pathways in glioblastoma and implications for combination therapy with temozolomide. *Neuro. Oncol.* **13**(4), 384-392 (2011).

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