

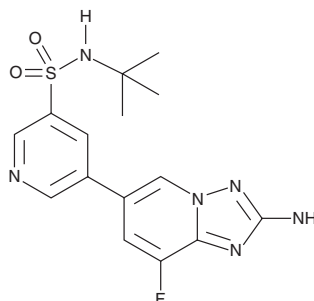
PRODUCT INFORMATION



CZC-24832

Item No. 14307

CAS Registry No.: 1159824-67-5
Formal Name: 5-(2-amino-8-fluoro[1,2,4]triazolo[1,5-a]pyridin-6-yl)-N-(1,1-dimethylethyl)-3-pyridinesulfonamide
MF: C₁₅H₁₇FN₆O₂S
FW: 364.4
Purity: ≥98%
UV/Vis.: λ_{max}: 218, 254, 292 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

CZC-24832 is supplied as a crystalline solid. A stock solution may be made by dissolving the CZC-24832 in the solvent of choice, which should be purged with an inert gas. CZC-24832 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of CZC-24832 in these solvents is approximately 5 mg/ml.

CZC-24832 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CZC-24832 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. CZC-24832 has a solubility of approximately 0.2 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

CZC-24832 is an inhibitor of PI3Kγ (IC₅₀ = 0.025 μM in a binding assay).¹ It is selective for PI3Kγ over PI3Kδ, PI3Kα, and PI3Kβ (IC₅₀s = 7.9, >10, and 1.26 μM, respectively, in binding assays), as well as a panel of 154 lipid and protein kinases.^{1,2} CZC-24832 inhibits phosphorylation of Akt induced by complement component 5a (C5a) in RAW 264.7 macrophages (IC₅₀ = 1.2 μM) and migration of isolated human granulocytes induced by fMLP (Item No. 21495; IC₅₀ = 1 μM).² It decreases disease severity in a mouse model of collagen-induced arthritis when administered at doses of 3 and 10 mg/kg.

References

1. Bell, K., Sunose, M., Ellard, K., *et al.* SAR studies around a series of triazolopyridines as potent and selective PI3Kγ inhibitors. *Bioorg. Med. Chem. Lett.* **22(16)**, 5257-5263 (2012).
2. Bergamini, G., Bell, K., Shimamura, S., *et al.* A selective inhibitor reveals PI3Kγ dependence of T(H)17 cell differentiation. *Nat. Chem. Biol.* **8(6)**, 576-582 (2012).

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, 11/07/2022

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