

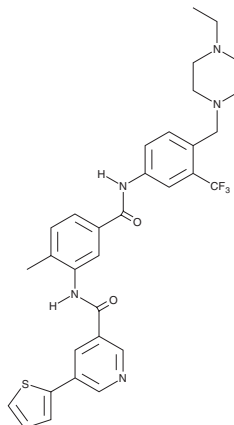
PRODUCT INFORMATION



ALW-II-41-27

Item No. 25275

CAS Registry No.: 1186206-79-0
Formal Name: N-[4-[(4-ethyl-1-piperazinyl)methyl]-3-(trifluoromethyl)phenyl]-4-methyl-3-[[[5-(2-thienyl)-3-pyridinyl]carbonyl]amino]-benzamide
MF: C₃₂H₃₂F₃N₅O₂S
FW: 607.7
Purity: ≥95%
UV/Vis.: λ_{max}: 277 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

ALW-II-41-27 is supplied as a crystalline solid. A stock solution may be made by dissolving the ALW-II-41-27 in the solvent of choice, which should be purged with an inert gas. ALW-II-41-27 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of ALW-II-41-27 in ethanol is approximately 0.5 mg/ml and approximately 10 mg/ml in DMSO and DMF.

ALW-II-41-27 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, ALW-II-41-27 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. ALW-II-41-27 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

ALW-II-41-27 is a multi-kinase inhibitor that inhibits EphB2, EphA3, Kit, FMS, VEGFR2/KDR, FLT1, FGR, Src, Lyn, BMX, and Bcr-Abl in tyrosine kinase-transformed Ba/F3 cells (EC₅₀s = <500 nM).¹ ALW-II-41-27 also inhibits DDR2 and Src (IC₅₀s = 51 and 14 nM, respectively) as well as wild-type and mutant RET kinases (IC₅₀s = 24.7, 94.2, and 15.8 nM for wild-type, RET^{V804L}, and RET^{V804M}, respectively).^{2,3} It reduces growth of NCI-H2286 and HCC-366 cancer cells (GI₅₀s = 0.51 and 0.65 μM, respectively) and RAT1 cells transformed by RET^{C634R} or RET^{M918T} (IC₅₀s = 44 and 56 nM, respectively). ALW-II-41-27 also inhibits growth of MDA-MB-231 breast cancer cells in a concentration-dependent manner and inhibits tumor growth *in vivo* in a mouse patient-derived xenograft (PDX) model of EphA2-overexpressing triple-negative breast cancer (TNBC).⁴

References

1. Choi, Y., Syeda, F., Walker, J.R., *et al.* Discovery and structural analysis of Eph receptor tyrosine kinase inhibitors. *Bioorg. Med. Chem. Lett.* **19(15)**, 4467-4470 (2009).
2. Terai, H., Tan, L., Beauchamp, E.M., *et al.* Characterization of DDR2 inhibitors for the treatment of DDR2 mutated nonsmall cell lung cancer. *ACS Chem. Biol.* **10(12)**, 2687-2696 (2015).
3. Moccia, M., Liu, Q., Guida, T., *et al.* Identification of novel small molecule inhibitors of oncogenic RET kinase. *PLoS One* **10(6)**, e0128364 (2015).
4. Song, W., Hwang, Y., Youngblood, V.M., *et al.* Targeting EphA2 impairs cell cycle progression and growth of basal-like/triple-negative breast cancers. *Oncogene* **36(40)**, 5620-5630 (2017).

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