

PRODUCT INFORMATION



AD57 (hydrochloride)

Item No. 13975

CAS Registry No.: 2320261-72-9
Formal Name: N-[4-[4-amino-1-(1-methylethyl)-1H-pyrazolo[3,4-d]pyrimidin-3-yl]phenyl]-N'-[3-(trifluoromethyl)phenyl]-urea, monohydrochloride

MF: C₂₂H₂₀F₃N₇O • HCl

FW: 491.9

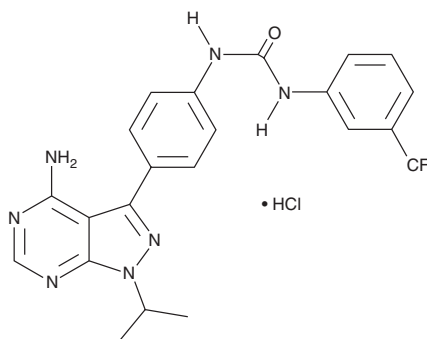
Purity: ≥98%

UV/Vis.: λ_{max}: 203, 271 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

AD57 (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the AD57 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. AD57 (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of AD57 (hydrochloride) in these solvents is approximately 5, 10, and 14 mg/ml, respectively.

AD57 (hydrochloride) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, AD57 (hydrochloride) should first be dissolved in DMF and then diluted with the aqueous buffer of choice. AD57 (hydrochloride) has a solubility of approximately 0.2 mg/ml in a 1:4 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Complex diseases, like cancer, may best be countered by compounds that affect multiple targets. AD57 is a polypharmacological cancer therapeutic, in that it is designed to modulate multiple targets related to cancer.¹ In *Drosophila*, it potently inhibits the receptor tyrosine kinase RET (IC₅₀ = 2 nM) and reduces the activity of numerous other kinases by more than 80% when given at 1 μM.¹ Most notably, AD57 interferes with kinases downstream of RET, including Src, Raf, and S6K, providing further efficacy in preventing signaling leading to invasion, proliferation, and metabolism relevant to cancer.¹

Reference

1. Dar, A.C., Das, T.K., Shokat, K.M., *et al.* Chemical genetic discovery of targets and anti-targets for cancer polypharmacology. *Nature* **486(7401)**, 80-84 (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, 04/01/2024

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