

Product Information

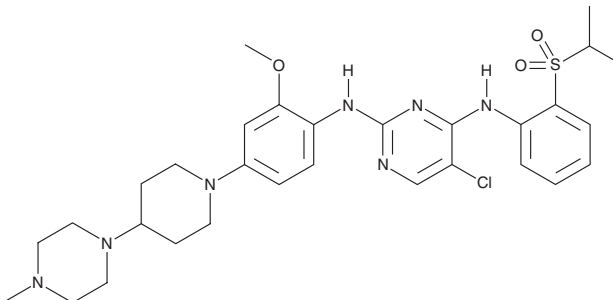


TAE684

Item No. 17670

CAS Registry No.: 761439-42-3
Formal Name: 5-chloro-N2-[2-methoxy-4-[4-(4-methyl-1-piperazinyl)-1-piperidinyl]phenyl]-N4-[2-[(1-methylethyl)sulfonyl]phenyl]-2,4-pyrimidinediamine

Synonym: NVP-TAE684
MF: C₃₀H₄₀ClN₇O₃S
FW: 614.2
Purity: ≥95%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid
UV/Vis.: λ_{max}: 209, 286 nm



Laboratory Procedures

For long term storage, we suggest that TAE684 be stored as supplied at -20°C. It should be stable for at least two years.

TAE684 is supplied as a crystalline solid. A stock solution may be made by dissolving the TAE684 in the solvent of choice. TAE684 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of TAE684 in these solvents is approximately 10 and 20 mg/ml, respectively.

TAE684 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, TAE684 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. TAE684 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Human anaplastic lymphoma kinase (ALK) is an oncogene that is amplified in neuroblastomas and when juxtaposed with various fusion partners, its constitutive kinase activity is associated with the development of a type of anaplastic large cell lymphoma (ALCL).^{1,2} TAE684 is an ALK inhibitor that blocks the proliferation of ALCL-derived and ALK-dependent cell lines with IC₅₀ values of 2-5 nM.¹ When tested against a panel of 35 cells transformed by various tyrosine kinases, TAE684 demonstrated 100- to 1,000-fold selectivity for inhibiting ALK-driven cell proliferation.¹ TAE684 treatment induces cell cycle arrest and apoptosis in ALK-dependent cell lines and has been used to suppress tumor growth in *in vivo* models of ALK-positive ALCL and neuroblastoma.^{1,2} TAE684 is also reported to inhibit the activity of the Parkinson's disease-linked leucine-rich repeat kinase 2 (IC₅₀s = 7.8 and 6.1 nM for wild-type and G2019S mutant LRRK2, respectively).³

References

1. Galkin, A.V., Melnick, J.S., Kim, S., *et al.* Identification of NVP-TAE684, a potent, selective, and efficacious inhibitor of NPM-ALK. *Proc. Natl. Acad. Sci. USA* **104**(1), 270-275 (2007).
2. Hasan, M.K., Nafady, A., Takatori, A., *et al.* ALK is a MYCN target gene and regulates cell migration and invasion in neuroblastoma. *Sci. Rep.* **3**, (2013).
3. Zhang, J., Deng, X., Choi, H.G., *et al.* Characterization of TAE684 as a potent LRRK2 kinase inhibitor. *Bioorg. Med. Chem. Lett.* **22**(55), 1864-1869 (2012).

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. 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

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery.**

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our Warranty and Limitation of Remedy located on our website and in our catalog.

Copyright Cayman Chemical Company, 06/30/2015

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com