

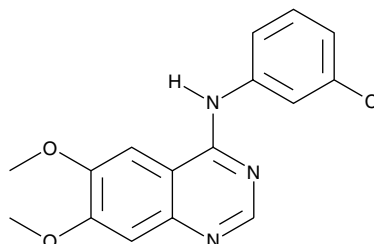
Product Information



AG-1478

Item No. 10010244

CAS Registry No.: 153436-53-4
Formal Name: N-(3-chlorophenyl)-6,7-dimethoxy-4-quinazolinamine
Synonyms: NSC 693255, Tyrphostin AG-1478
MF: C₁₆H₁₄ClN₃O₂
FW: 315.8
Purity: ≥98%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid
UV/Vis.: λ_{max}: 222, 255, 346 nm



Laboratory Procedures

For long term storage, we suggest that AG-1478 be stored as supplied at -20°C. It should be stable for at least two years.

AG-1478 is supplied as a crystalline solid. A stock solution may be made by dissolving the AG-1478 in an organic solvent purged with an inert gas. AG-1478 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of AG-1478 in ethanol, DMSO, and DMF is approximately 0.5, 1.5, and 1 mg/ml, respectively.

AG-1478 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, AG-1478 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. AG-1478 has a solubility of approximately 0.1 mg/ml in a 1:8 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Protein tyrosine kinase (PTK) inhibitors are potential antiproliferative agents for diseases caused by the hyperactivity of PTK. Tyrphostins are a class of antiproliferative compounds which act as PTK blockers. PTK inhibitors which preferentially inhibit the epidermal growth factor receptor (EGFR) kinase and block EGFR-dependent cell proliferation. AG-1478 is an inhibitor of EGFR kinase with an IC₅₀ value of 3 nM.¹ Due to its potency and selectivity, AG-1478 has been used in a broad range of studies. It reversibly inhibits rat brain Kv1.5 potassium channels (IC₅₀ = 9.8 μM) independent of PTK activity.² AG-1478 also inhibits the growth of leiomyoma and myometrium cell cultures with IC₅₀ values of 5.6 and 5.7 μM, respectively.³ This inhibitor suppresses MAP kinase activation and strongly inhibits induction of *fos* gene expression and DNA synthesis.⁴

References

- Levitzi, A. and Gazit, A. Tyrosine kinase inhibition: An approach to drug development. *Science* **267**(5205), 1782-1788 (1995).
- Choi, B.H., Choi, J.-S., Rhie, D.-J., *et al.* Direct inhibition of the cloned Kv1.5 channel by AG-1478, a tyrosine kinase inhibitor. *Am. J. Physiol. Cell Physiol.* **282**, c1461-c1468 (2002).
- Shushan, A., Rojansky, N., Laufer, N., *et al.* The AG1478 tyrosine kinase inhibitor is an effective suppressor of leiomyoma cell growth. *Hum. Reprod.* **19**(9), 1957-1967 (2004).
- Daub, H., Weiss, F.U., Wallasch, C., *et al.* Role of transactivation of the EGF receptor in signalling by G-protein-coupled receptors. *Nature* **379**, 557-560 (1996).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/10010244

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, 04/15/2015

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