

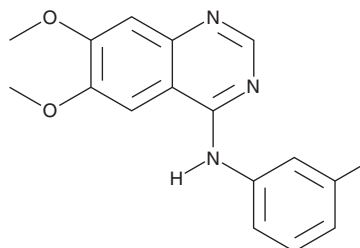
PRODUCT INFORMATION



AG-1557

Item No. 17989

CAS Registry No.: 189290-58-2
Formal Name: N-(3-iodophenyl)-6,7-dimethoxy-4-quinazolinamine
Synonym: Tyrphostin AG-1557
MF: C₁₆H₁₄IN₃O₂
FW: 407.2
Purity: ≥95%
UV/Vis.: λ_{max}: 221, 252, 345 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

AG-1557 is supplied as a crystalline solid. A stock solution may be made by dissolving the AG-1557 in the solvent of choice, which should be purged with an inert gas. AG-1557 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of AG-1557 in these solvents is approximately 0.5, 1.5, and 1 mg/ml, respectively.

AG-1557 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, AG-1557 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. AG-1557 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.

Description

AG-1557 is an inhibitor of epidermal growth factor receptor (EGFR) tyrosine kinase with a pIC₅₀ value of 8.194.¹

Reference

1. Joshi, A., Gadhwal, M., and Joshi, U.J. Identification of potential novel EGFR inhibitors using a combination of pharmacophore and docking methods. *Int. J. Pharm. Pharmaceut. Sci.* **7(6)** (2015).

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, 11/08/2022

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