

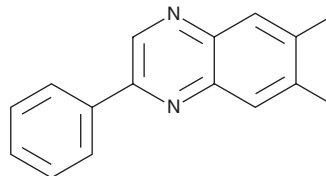
PRODUCT INFORMATION



AG-1295

Item No. 14529

CAS Registry No.: 71897-07-9
Formal Name: 6,7-dimethyl-2-phenyl-quinoxaline
Synonyms: NSC 380341, Tyrphostin AG-1295
MF: C₁₆H₁₄N₂
FW: 234.3
Purity: ≥98%
UV/Vis.: λ_{max}: 263, 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-1295 is supplied as a crystalline solid. A stock solution may be made by dissolving the AG-1295 in the solvent of choice, which should be purged with an inert gas. AG-1295 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of AG-1295 in these solvents is approximately 2.5, 2, and 5 mg/ml, respectively.

AG-1295 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, AG-1295 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. AG-1295 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

AG-1295 is a quinoxaline-type tyrphostin that acts as a potent and selective inhibitor of PDGF receptor kinase *in vitro* and in Swiss 3T3 cells (IC₅₀s range from 0.3-1 μM).¹ It inhibits PDGF-stimulated DNA synthesis with an IC₅₀ value of 2.5 μM without affecting activity of the EGF receptor.¹

Reference

1. Kovalenko, M., Gazit, A., Böhmer, A., *et al.* Selective platelet-derived growth factor receptor kinase blockers reverse sis-transformation. *Cancer Res.* **54(23)**, 6106-6114 (1994).

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, 12/15/2022

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