

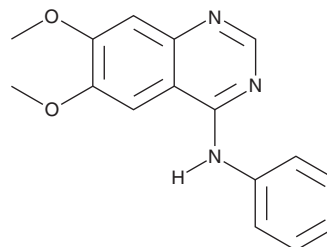
PRODUCT INFORMATION



WHI-P258

Item No. 23453

CAS Registry No.: 21561-09-1
Formal Name: 6,7-dimethoxy-N-phenyl-4-quinazolinamine
MF: C₁₆H₁₅N₃O₂
FW: 281.3
Purity: ≥98%
UV/Vis.: λ_{max}: 201, 223, 254, 346 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

WHI-P258 is supplied as a crystalline solid. A stock solution may be made by dissolving the WHI-P258 in the solvent of choice, which should be purged with an inert gas. WHI-P258 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of WHI-P258 in these solvents is approximately 15 and 25 mg/ml, respectively.

WHI-P258 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, WHI-P258 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. WHI-P258 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

WHI-P258 is a quinazoline compound that modeling studies suggested would bind to the active site of JAK3 with an estimated K_i value of 72 μM.¹ However, it is inactive at JAK3 (IC₅₀ = >300 μM) and has been used as a negative control for structurally similar compounds that inhibit platelet aggregation and herpes simplex virus 1 (HSV-1) replication.^{2,3}

References

1. Sudbeck, E.A., Liu, X.P., Narla, R.K., *et al.* Structure-based design of specific inhibitors of janus kinase 3 as apoptosis-inducing antileukemic agents. *Clin. Cancer Res.* **5(6)**, 1569-1582 (1999).
2. Yakota, S.-i., Yokosawa, N., Okabayashi, T., *et al.* Induction of suppressor of cytokine signaling-3 by herpes simplex virus type 1 confers efficient viral replication. *Virology* **338(1)**, 173-181 (2005).
3. Tibbles, H.E., Vassilev, A., Wendorf, H., *et al.* Role of a JAK3-dependent biochemical signaling pathway in platelet activation and aggregation. *J. Biol. Chem.* **276(21)**, 17815-17822 (2001).

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, 11/15/2022

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