

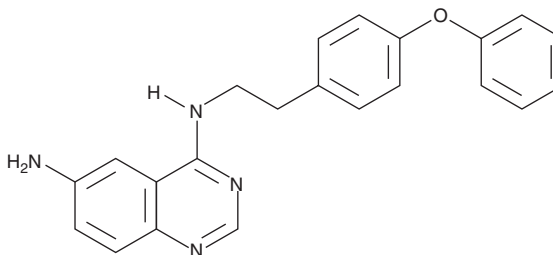
PRODUCT INFORMATION



QNZ

Item No. 10006734

CAS Registry No.: 545380-34-5
Formal Name: N4-[2-(4-phenoxyphenyl)ethyl]-4,6-quinazolinediamine
Synonym: CAY10470
MF: C₂₂H₂₀N₄O
FW: 356.4
Purity: ≥98%
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

QNZ is supplied as a crystalline solid. A stock solution may be made by dissolving the QNZ in the solvent of choice, which should be purged with an inert gas. QNZ is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of QNZ in these solvents is approximately 10, 20, and 30 mg/ml, respectively.

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

Nuclear factor- κ B (NF- κ B) is an ubiquitous transcription factor that functions to enhance the transcription of pro-inflammatory cytokines like TNF α , Interleukin-1 β , IL-6, and IL-8.¹ QNZ is an inhibitor of NF- κ B activation with an IC₅₀ of 11 nM in human Jurkat cells. It also inhibits TNF- α production from lipopolysaccharide-stimulated mouse splenocytes with an IC₅₀ of 7 nM. QNZ showed efficacy *in vivo* by reducing edema formation in carrageenin-induced inflammation of rat hind paw.²

References

1. Karin, M., Yamamoto, Y., and Wang, Q.M. The IKK NF- κ B system: A treasure trove for drug development. *Nat. Rev. Drug Discov.* **3**(1), 17-26 (2004).
2. Tobe, M., Isobe, Y., Tomizawa, H., *et al.* Discovery of quinazolines as a novel structural class of potent inhibitors of NF- κ B activation. *Bioorg. Med. Chem.* **11**(3), 383-391 (2003).

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, 11/22/2022

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