

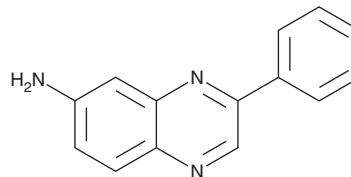
PRODUCT INFORMATION



PAQ

Item No. 20234

CAS Registry No.: 943902-10-1
Formal Name: 3-phenyl-6-quinoxalinamine
Synonym: 3-Phenyl-6-aminoquinoxaline
MF: C₁₄H₁₁N₃
FW: 221.3
Purity: ≥95%
UV/Vis.: λ_{max}: 248, 282, 403 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

PAQ is supplied as a crystalline solid. A stock solution may be made by dissolving the PAQ in the solvent of choice, which should be purged with an inert gas. PAQ is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of PAQ in ethanol is approximately 2 mg/ml and approximately 10 mg/ml in DMSO and DMF.

PAQ is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, PAQ should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. PAQ has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

PAQ is a neuroprotective agent.¹ It protects dopaminergic neurons from cell death without inhibiting glial cell proliferation in a rat midbrain culture model of Parkinson's disease when used at a concentration of 10 μM. PAQ (25 or 50 mg/kg, twice per day) prevents loss of dopaminergic neurons in the substantia nigra in a mouse model of Parkinson's disease induced by MPTP.

Reference

1. Le Douaron, G., Ferrié, L., Sepulveda-Diaz, J.E., *et al.* New 6-aminoquinoxaline derivatives with neuroprotective effect on dopaminergic neurons in cellular and animal parkinson disease models. *J. Med. Chem.* **59**(13), 6169-6186 (2016).

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, 02/06/2025

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