

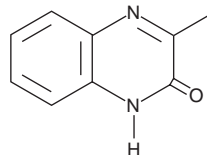
PRODUCT INFORMATION



3-methyl-2-Quinoxalinone

Item No. 13006

CAS Registry No.: 14003-34-0
Formal Name: 3-methyl-2(1H)-quinoxalinone
Synonyms: NSC 17798, NSC 38589, NSC 149824
MF: C₉H₈N₂O
FW: 160.1
Purity: ≥98%
UV/Vis.: λ_{max}: 203, 229, 278, 339 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

3-methyl-2-Quinoxalinone is supplied as a crystalline solid. A stock solution may be made by dissolving the 3-methyl-2-quinoxalinone in the organic solvent of choice, which should be purged with an inert gas. 3-methyl-2-Quinoxalinone is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of 3-methyl-2-quinoxalinone in ethanol is approximately 1 mg/ml and approximately 20 mg/ml in DMSO and DMF.

3-methyl-2-Quinoxalinone is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 3-methyl-2-quinoxalinone should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. 3-methyl-2-Quinoxalinone 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

3-methyl-2-Quinoxalinone is a synthetic intermediate useful for pharmaceutical synthesis.

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, 12/13/2022

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