

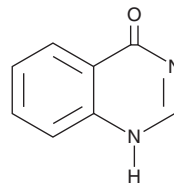
PRODUCT INFORMATION



4-HQN

Item No. 32821

CAS Registry No.: 491-36-1
Formal Name: 4(3H)-quinazolinone
Synonym: 4-Hydroxyquinazoline
MF: $C_8H_6N_2O$
FW: 146.1
Purity: $\geq 97\%$
UV/Vis.: λ_{\max} : 223 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

4-HQN is supplied as a crystalline solid. A stock solution may be made by dissolving the 4-HQN in the solvent of choice, which should be purged with an inert gas. 4-HQN is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of 4-HQN in these solvents is approximately 1 mg/ml.

4-HQN is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 4-HQN should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. 4-HQN has a solubility of approximately 0.09 mg/ml in a 1:10 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

4-HQN is a heterocyclic building block.^{1,2} It has been used in the synthesis of compounds with antimalarial or antimicrobial activities *in vitro*.

References

1. Kikuchi, H., Yamamoto, K., Horoiwa, S., *et al.* Exploration of a new type of antimalarial compounds based on febrifugine. *J. Med. Chem.* **49**(15), 4698-4706 (2006).
2. Yan, B.-R., Lv, X.-Y., Du, H., *et al.* Synthesis and biological activities of novel quinazolinone derivatives containing a 1,2,4-triazolylthioether moiety. *De Gruyter* **70**(7), 983-993 (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, 11/10/2022

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