

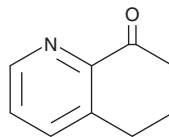
PRODUCT INFORMATION



6,7-Dihydro-5H-quinolin-8-one

Item No. 30480

CAS Registry No.: 56826-69-8
Formal Name: 6,7-dihydro-8(5H)-quinolinone
Synonyms: 8-Aza-1-tetralone, 5,6-Dihydro-8(7H)-quinolinone
MF: C₉H₉NO
FW: 147.2
Purity: ≥98%
UV/Vis.: λ_{max}: 230, 282 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

6,7-Dihydro-5H-quinolin-8-one is supplied as a crystalline solid. A stock solution may be made by dissolving the 6,7-dihydro-5H-quinolin-8-one in the solvent of choice, which should be purged with an inert gas. 6,7-Dihydro-5H-quinolin-8-one is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of 6,7-dihydro-5H-quinolin-8-one in these solvents is approximately 30 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of 6,7-dihydro-5H-quinolin-8-one can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of 6,7-dihydro-5H-quinolin-8-one in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

6,7-Dihydro-5H-quinolin-8-one is a synthetic intermediate.^{1,2} It has been used in the synthesis of tetrahydropyridoazepinones and thiosemicarbazones with anticancer activity.

References

1. Lemke, T.L., Shek, T.W., Cates, L.A., *et al.* Synthesis of 5,6-dihydro-8(7H)-quinolinone thiosemicarbazones as potential antitumor agents. *J. Med. Chem.* **20(10)**, 1351-1354 (1977).
2. Jössang-Yanagida, A. and Gansser, C. Tetrahydropyridoazepines and tetrahydropyridoazepinones from the corresponding dihydroquinolones. *J. Heterocycl. Chem.* **15(2)**, 249-251 (1977).

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, 10/27/2022

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