

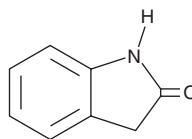
PRODUCT INFORMATION



Oxindole

Item No. 23851

CAS Registry No.: 59-48-3
Formal Name: 1,3-dihydro-2H-indol-2-one
Synonyms: 2-Indolone, NSC 274863
MF: C₈H₇NO
FW: 133.1
Purity: ≥98%
UV/Vis.: λ_{max}: 249 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

Oxindole is supplied as a crystalline solid. A stock solution may be made by dissolving the oxindole in the solvent of choice. Oxindole is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of oxindole in ethanol and DMF is approximately 10 mg/ml and approximately 3 mg/ml in DMSO.

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

Description

Oxindole is an aromatic heterocycle and starting material in the synthesis of numerous biologically active compounds including, but not limited to, anticancer, antibacterial, and antidiabetic agents, enzyme inhibitors, and channel blockers.^{1,2}

References

1. Ziarani, G.M., Gholamzadeh, P., Lashgari, N., *et al.* Oxindole as starting material in organic synthesis. *ARKIVOC* **2013**(1), 470-535 (2013).
2. Kaur, M., Singh, M., Chadha, N., *et al.* Oxindole: A chemical prism carrying plethora of therapeutic benefits. *Eur. J. Med. Chem.* **123**, 858-894 (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, 12/22/2022

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