

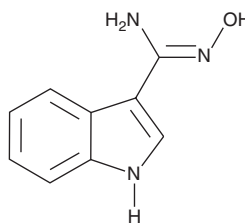
PRODUCT INFORMATION



Indole-3-amidoxime

Item No. 11060

CAS Registry No.: 95649-37-9
Formal Name: N-hydroxy-1H-indole-3-carboximidamide
MF: $C_9H_9N_3O$
FW: 175.2
Purity: $\geq 97\%$
UV/Vis.: λ_{max} : 216, 280 nm
Supplied as: A crystalline solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Indole-3-amidoxime is supplied as a crystalline solid. A stock solution may be made by dissolving the indole-3-amidoxime in an organic solvent purged with an inert gas. Indole-3-amidoxime is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of indole-3-amidoxime in these solvents is approximately 14 and 12.5 mg/ml, respectively.

Indole-3-amidoxime is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, indole-3-amidoxime should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Indole-3-amidoxime has a solubility of approximately 0.1 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

Indole-3-amidoxime 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, 12/12/2022

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