

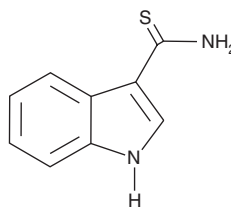
# PRODUCT INFORMATION



## Indole-3-thio Carboxamide

Item No. 11065

**CAS Registry No.:** 59108-90-6  
**Formal Name:** 1H-indole-3-carbothioamide  
**MF:** C<sub>9</sub>H<sub>8</sub>N<sub>2</sub>S  
**FW:** 176.2  
**Purity:** ≥97%  
**UV/Vis.:** λ<sub>max</sub>: 213, 252, 276, 327 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

Indole-3-thio Carboxamide is supplied as a crystalline solid. A stock solution may be made by dissolving the indole-3-thio carboxamide in the solvent of choice, which should be purged with an inert gas. Indole-3-thio Carboxamide is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of indole-3-thio carboxamide in ethanol and DMSO is approximately 20 mg/ml and approximately 30 mg/ml in DMF.

Indole-3-thio Carboxamide is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, indole-3-thio carboxamide should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Indole-3-thio Carboxamide has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Indole-3-thio Carboxamide 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