

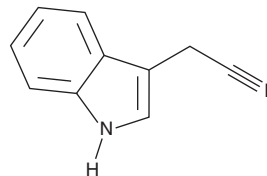
PRODUCT INFORMATION



Indole-3-acetonitrile

Item No. 38751

CAS Registry No.: 771-51-7
Formal Name: 1H-indole-3-acetonitrile
Synonyms: IAN, 3-Indoleacetonitrile, 3-Indolylacetonitrile, NSC 523272
MF: C₁₀H₈N₂
FW: 156.2
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Synthetic



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

Laboratory Procedures

Indole-3-acetonitrile is supplied as a solid. A stock solution may be made by dissolving the indole-3-acetonitrile in the solvent of choice, which should be purged with an inert gas. Indole-3-acetonitrile is soluble in methanol.

Description

Indole-3-acetonitrile is an indole alkaloid that has been found in cabbage and has antibacterial activity.^{1,2} It reduces biofilm formation by *E. coli* and *P. aeruginosa* when used at a concentration of 100 µg/ml.¹ The levels of indole-3-acetonitrile are increased in *P. brassicae*-infected cabbage roots.²

References

1. Lee, J.-H., Cho, M.H., and Lee, J. 3-indolylacetonitrile decreases *Escherichia coli* O157:H7 biofilm formation and *Pseudomonas aeruginosa* virulence. *Environ. Microbiol.* **13(1)**, 62-73 (2011).
2. Kavanagh, J.A. and Williams, P.H. Indole auxins in *Plasmodiophora* infected cabbage roots and hypocotyls. *Trans. Br. Mycol. Soc.* **77(1)**, 125-129 (1981).

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, 12/06/2023

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