

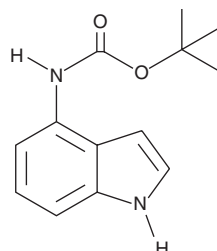
PRODUCT INFORMATION



(1H-Indol-4-yl)-Carbamic Acid *tert*-butyl ester

Item No. 10007087

CAS Registry No.: 819850-13-0
Formal Name: 1H-indol-4-yl-carbamic acid, 1,1-dimethylethyl ester
MF: C₁₃H₁₆N₂O₂
FW: 232.3
Purity: ≥95%
UV/Vis.: λ_{max}: 219, 278 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

(1H-Indol-4-yl)-Carbamic acid *tert*-butyl ester is supplied as a crystalline solid. A stock solution may be made by dissolving the (1H-Indol-4-yl)-carbamic acid *tert*-butyl ester in the solvent of choice, which should be purged with an inert gas. (1H-Indol-4-yl)-Carbamic acid *tert*-butyl ester is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of (1H-Indol-4-yl)-carbamic acid *tert*-butyl ester in these solvents is approximately 5, 25, and 30, respectively.

(1H-Indol-4-yl)-Carbamic acid *tert*-butyl ester is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, (1H-Indol-4-yl)-carbamic acid *tert*-butyl ester should first be dissolved in DMF and then diluted with the aqueous buffer of choice. (1H-Indol-4-yl)-Carbamic acid *tert*-butyl ester has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

(1H-Indol-4-yl)-Carbamic Acid *tert*-butyl ester 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, 12/10/2022

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