

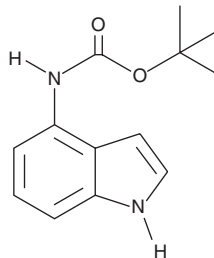
# Product Information



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

Catalog No. 10007087

**CAS Registry No.:** 819850-13-0  
**Formal Name:** 1H-indol-4-yl-carbamic acid, 1,1-dimethylethyl ester  
**MF:** C<sub>13</sub>H<sub>16</sub>N<sub>2</sub>O<sub>2</sub>  
**FW:** 232.3  
**Purity:** ≥95%  
**Stability:** ≥2 years at -20°C  
**Supplied as:** A crystalline solid



### Applications

For long term storage, we suggest that (1H-Indol-4-yl)-carbamic acid *tert*-butyl ester be stored as supplied at -20°C. It should be stable for at least two years.

(1H-Indol-4-yl)-Carbamic acid *tert*-butyl ester is a synthetic intermediate useful for pharmaceutical synthesis.

### Related Products

4-(1-piperazinyl)-1H-Indole (dihydrochloride) - Cat. No. 10007075 • CAY10491 - Cat. No. 10007076 • CAY10492 - Cat. No. 10007077 • 4-Cyanoindole - Cat. No. 10007084 • 4-Aminomethylindole - Cat. No. 10007085 • 4-formyl Indole - Cat. No. 10007086 • CAY10494 - Cat. No. 10007126 • 7-Piperazin-1-yl-thieno[2,3-c] Pyridine (hydrochloride) - Cat. No. 10007136 • 8-Piperazin-1-yl-Isoquinoline (hydrochloride) - Cat. No. 10007146 • 7-piperazin-1-yl-Isoquinoline (hydrochloride) - Cat. No. 10007147 • 6-piperazin-1-yl-Isoquinoline (hydrochloride) - Cat. No. 10007148

### Cayman Chemical

**Mailing address**  
1180 E. Ellsworth Road  
Ann Arbor, MI  
48108 USA

**Phone**  
(800) 364-9897  
(734) 971-3335

**Fax**  
(734) 971-3640

**E-Mail**  
custserv@caymanchem.com

**Web**  
www.caymanchem.com

**WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.**

#### MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent *via* email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will meet our specifications at the time of delivery.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a refund of the purchase price, or at Cayman's option, the replacement, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy** located on our website and in our catalog.

Copyright Cayman Chemical Company, 03/04/2010