

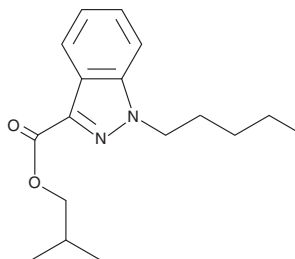
Product Information



Isobutyl 1-pentyl-1H-indazole-3-carboxylate RM

Item No. 17522

Formal Name: isobutyl 1-pentyl-1H-indazole-3-carboxylate
MF: C₁₇H₂₄N₂O₂
FW: 288.4
Purity: ≥98%
Storage: -20°C
Supplied as: A neat oil



Description

NNEI 2'-indazole isomer (Item No. 17520) is an analog of the synthetic cannabinoid NNEI (Item No. 15001) that has the alkyl group at the two position of an indazole scaffold.¹ Isobutyl 1-pentyl-1H-indazole-3-carboxylate is a compound recently identified in cannabinoid mixtures in Japan.¹ It is considered to be a by-product of the preparation of NNEI 2'-indazole isomer.¹ The physiological and toxicological properties of this compound are not known. This product is intended for forensic and research applications.

Cayman Chemical products containing Reference Material (RM) in the name are highly qualified materials manufactured and tested to meet ISO/IEC 17025:2005 and ISO Guide 34:2009 international guidelines for reference materials. These materials are tested using validated analytical methods on qualified instrumentation to ensure traceability to SI units of measure.

Reference

1. Uchiyama, N., Shimokawa, Y., Kikura-Hanajiri, R., *et al.* A synthetic cannabinoid FDU-NNEI, two 2H-indazole isomers of synthetic cannabinoids AB-CHMINACA and NNEI indazole analog (MN-18), a phenethylamine derivative N-OH-EDMA, and a cathinone derivative dimethoxy- α -PHP, newly identified in illegal products. *Forensic Toxicol.* (2015).

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.

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 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, 06/29/2015