

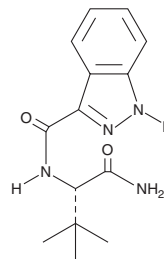
PRODUCT INFORMATION



ADB-INACA

Item No. 37729

CAS Registry No.: 1887742-42-8
Formal Name: N-[(1S)-1-(aminocarbonyl)-2,2-dimethylpropyl]-1H-indazole-3-carboxamide
MF: C₁₄H₁₈N₄O₂
FW: 274.3
Purity: ≥98%
UV/Vis.: λ_{max}: 295 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥5 years



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

Description

AADB-INACA (Item No. 37729) is an analytical reference standard categorized as a synthetic cannabinoid metabolite and precursor in the synthesis of synthetic cannabinoids.¹ ADB-INACA is a metabolite of ADB-PINACA (Item Nos. ISO60212 | ISO00150) and 5-fluoro ADB-PINACA (Item No. 14764) and a precursor in the synthesis of synthetic cannabinoids, including ADB-FUBINACA (Item No. 14292) and MAB-CHMINACA (Item No. 16616). This product is intended for research and forensic applications.

Reference

1. Carlier, J., Diao, X., Scheidweiler, K.B., *et al.* Distinguishing intake of new synthetic cannabinoids ADB-PINACA and 5F-ADB-PINACA with human hepatocyte metabolites and high-resolution mass spectrometry. *Clin. Chem.* **63**(5), 1008-1021 (2017).

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, 02/27/2024

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