PRODUCT INFORMATION



AB-CHMINACA 2'-indazole isomer

Item No. 17521

CAS Registry No.:	2237296-99-8	
Formal Name:	N-[(1S)-1-(aminocarbonyl)-2-	
	methylpropyl]-2-(cyclohexylmethyl)-2H-	
	indazole-3-carboxamide	
MF:	C ₂₀ H ₂₈ N ₄ O ₂	O. N
FW:	356.5	
Purity:	≥98%	Ń,
Supplied as:	A neat solid	H ₂ N T T
Storage:	-20°C	
Stability:	≥2 years	
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysi		

rovided on each certificate of analysis.

Description

AB-CHMINACA (Item No. 15434) is a synthetic cannabinoid (CB) that is structurally related to AB-FUBINACA (Item No. 14039), a potent agonist of the central CB₁ receptor.^{1,2} AB-CHMINACA 2'-indazole isomer is an isomer of AB-CHMINACA in which the cyclohexyImethyl moiety is attached at the 2 position of the indazole group, rather than at the 1 position.³ The physiological and toxicological properties of this compound are not known. This product is intended for forensic and research applications.

This product is qualified as a Reference Material that has been manufactured and tested to ISO/IEC 17025 and ISO 17034 international standards.

References

- 1. Uchiyama, N., Matsuda, S., Wakana, D., et al. New cannabimimetic indazole derivatives, N-(1-amino-3methyl-1-oxobutan-2-yl)-1-pentyl-1H-indazole-3-carboxamide (AB-PINACA) and N-(1-amino-3methyl-1-oxobutan-2-yl)-1-(4-fluorobenzyl)-1H-indazole-3-carboxamide (AB-FUBINACA) identified as designer drugs in illegal products. Forensic Toxicol. 31(1), 93-100 (2013).
- 2. Buchler, I.P., Hayes, M.J., Hedge, S.G., et al. Indazole derivatives. WO 2009/106982 A1 (2009), 1-283, PCT/IB2009/000432.
- 3. Uchiyama, N., Shimokawa, Y., Kikura-Hanajiri, R., et al. Asynthetic 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. 33(2), 244-259 (2015).

WARNING THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFFTY 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.

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, 08/04/2022

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