

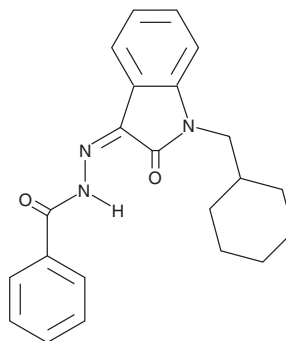
PRODUCT INFORMATION



BZO-CHMOXIZID

Item No. 34956

CAS Registry No.: 1048973-67-6
Formal Name: benzoic acid, (2Z)-2-[1-(cyclohexylmethyl)-1,2-dihydro-2-oxo-3H-indol-3-ylidene]hydrazide
Synonym: CHM-MDA-19
MF: C₂₂H₂₃N₃O₂
FW: 361.4
Purity: ≥98%
UV/Vis.: λ_{max}: 203, 277, 343 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥5 years
Item Origin: Synthetic



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

Description

BZO-CHMOXIZID (Item No. 34956) is an analytical reference standard categorized as a synthetic cannabinoid.^{1,2} This product is intended for research and forensic applications.

References

1. Deventer, M.H., Van Uytvanghe, K., Vinckier, I.M.J., *et al.* Cannabinoid receptor activation potential of the next generation, generic ban evading OXIZID synthetic cannabinoid receptor agonists. *Drug Test Anal.* (2022).
2. Diaz, P., Xu, J., Astruc-Diaz, F., *et al.* Design and synthesis of a novel series of N-alkyl isatin acylhydrazone derivatives that act as selective cannabinoid receptor 2 agonists for the treatment of neuropathic pain. *J. Med. Chem.* **51**(16), 4932-4947 (2008).

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, 06/29/2023

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