

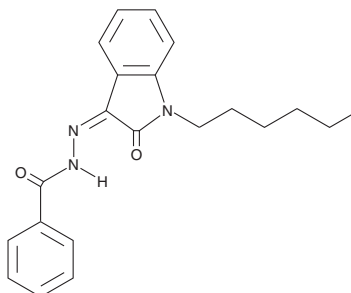
# PRODUCT INFORMATION



## MDA 19

Item No. 10563

**CAS Registry No.:** 1048973-47-2  
**Formal Name:** (2Z)-2-(1-hexyl-1,2-dihydro-2-oxo-3H-indol-3-ylidene)hydrazide, benzoic acid  
**Synonym:** BZO-HEXOXIZID  
**MF:** C<sub>21</sub>H<sub>23</sub>N<sub>3</sub>O<sub>2</sub>  
**FW:** 349.4  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 276, 342 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥5 years



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

### Description

MDA 19 (Item No. 10563) is an analytical reference standard categorized as a synthetic cannabinoid.<sup>1</sup> MDA 19 has been found in illicit substances seized by law enforcement. It reduces tactile allodynia in rats and mice.<sup>2,3</sup> 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).
3. Xu, J.J., Diaz, P., Astruc-Diaz, F., *et al.* Pharmacological characterization of a novel cannabinoid ligand, MDA19, for treatment of neuropathic pain. *Anesth. Analg.* **111(1)**, 99-109 (2010).

#### 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, 07/18/2023

#### 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](http://WWW.CAYMANCHEM.COM)