

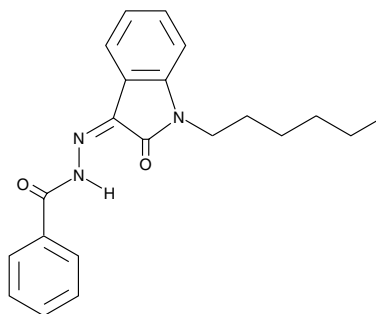
# Product Information



## MDA 19

Item No. 10563

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



### Laboratory Procedures

For long term storage, we suggest that MDA 19 be stored as supplied at -20°C. It should be stable for at least two years.

MDA 19 is supplied as a crystalline solid. A stock solution may be made by dissolving the MDA 19 in the solvent of choice. MDA 19 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of MDA 19 in these solvents is approximately 3, 10, and 20 mg/ml, respectively.

If aqueous stock solutions are required for biological experiments, they can best be prepared by diluting the organic solvent into aqueous buffers or isotonic saline. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

MDA 19 is a selective agonist of the human peripheral cannabinoid (CB<sub>2</sub>) receptor, with an EC<sub>50</sub> value for activation (63.4 nM) that is 14-fold lower than that for central cannabinoid (CB<sub>1</sub>) activation (EC<sub>50</sub> = 867 nM).<sup>1</sup> Surprisingly, it acts as an inverse agonist at the rat CB<sub>2</sub> receptor in cell-based functional assays.<sup>2</sup> MDA 19 attenuates tactile allodynia produced by spinal nerve ligation or paclitaxel in a dose-related manner in rats and in CB<sub>2</sub><sup>+/+</sup> mice but not in CB<sub>2</sub><sup>-/-</sup> mice, indicating that CB<sub>2</sub> receptors mediated the effects of MDA 19.<sup>1,2</sup> These effects of MDA 19 are blocked by the selective CB<sub>2</sub> antagonist AM630.<sup>1</sup>

### References

1. 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**, 4932-4947 (2008).
2. 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**, 99-109 (2010).

### Related Products

For a list of related products please visit: [www.caymanchem.com/catalog/10563](http://www.caymanchem.com/catalog/10563)

**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, 01/07/2014

### 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](mailto:custserv@caymanchem.com)

#### Web

[www.caymanchem.com](http://www.caymanchem.com)