

PRODUCT INFORMATION

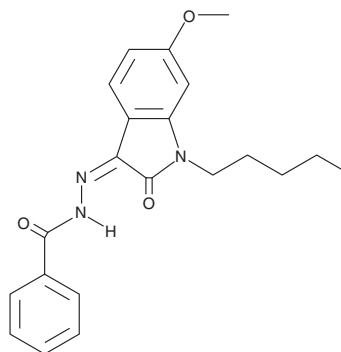


MDA 77

Item No. 10639

CAS Registry No.: 1103774-21-5
Formal Name: 2Z-(1,2-dihydro-6-methoxy-2-oxo-1-pentyl-3H-indol-3-ylidene)hydrazide benzoic acid

MF: C₂₁H₂₃N₃O₃
FW: 365.4
Purity: ≥98%
UV/Vis.: λ_{max}: 284, 353 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



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

Description

MDA 77 is a selective inverse agonist of the human peripheral cannabinoid (CB₂) receptor that demonstrates an EC₅₀ value of 5.8 nM at CB₂ and no activity at central cannabinoids (CB₁) receptor in functional binding assays.¹ The *in vivo* activity of this compound has not been reported. This product is intended for research and forensic applications.

Reference

1. Diaz, P., Phatuk, S.S., Xu, J., *et al.* 6-Methoxy-N-alkyl isatin acylhydrazone derivatives as a novel series of potent selective cannabinoid receptor 2 inverse agonists: Design, synthesis, and binding mode prediction. *J. Med. Chem.* **52**(2), 433-44 (2009).

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, 08/15/2022

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