

PRODUCT INFORMATION

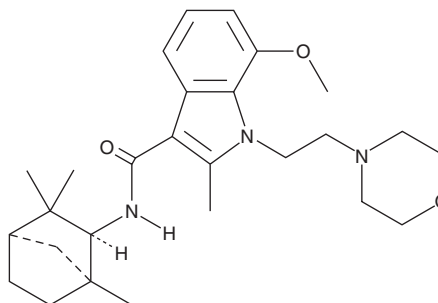


MN-25-2-methyl derivative

Item No. 9001455

CAS Registry No.: 501927-29-3
Formal Name: 7-methoxy-2-methyl-1-[2-(4-morpholinyl)ethyl]-N-[(1S,2S,4R)-1,3,3-trimethylbicyclo[2.2.1]hept-2-yl]-1H-indole-3-carboxamide

MF: C₂₇H₃₉N₃O₃
FW: 453.6
Purity: ≥95%
UV/Vis.: λ_{max}: 223, 284 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

MN-25-2-methyl derivative is a cannabinoid (CB) receptor ligand with K_i values of 8 and 29 nM at the central CB₁ receptor and peripheral CB₂ receptor, respectively.¹ This product is intended for research and forensic applications.

Reference

1. Wroblewski, S.T., Chen, P., Hynes, J., Jr., *et al.* Rational design and synthesis of an orally active indolopyridone as a novel conformationally constrained cannabinoid ligand possessing antiinflammatory properties. *J. Med. Chem.* **46**(11), 2110-2116 (2003).

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, 07/11/2023

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