

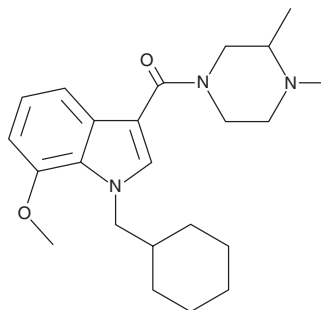
PRODUCT INFORMATION



(±)-ORG 28611

Item No. 15266

CAS Registry No.: 784138-08-5
Formal Name: [1-(cyclohexylmethyl)-7-methoxy-1H-indol-3-yl](3,4-dimethyl-1-piperazinyl)-methanone
MF: C₂₃H₃₃N₃O₂
FW: 383.5
Purity: ≥98%
UV/Vis.: λ_{max}: 219, 288 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

(±)-ORG 28611 is the racemic mixture of the water soluble, full agonist of the human cannabinoid (CB) receptors central CB₁ (pEC₅₀ = 7.6; pK_i = 8.9) and peripheral CB₂ (pK_i = 8.8).¹ ORG 28611 (0.1-1 μM/kg, intravenous) demonstrated antinociceptive activity in the mouse tail flick test and the mouse formalin paw test, and it was further evaluated clinically for use as a potential intravenous analgesic agent.¹ This product is intended for forensic and research applications.

References

1. Adam, J.M., Cairns, J., Caulfield, W., *et al.* Design, synthesis, and structure-activity relationships of indole-3-carboxamides as novel water soluble cannabinoid CB₁ receptor agonists. *Med. Chem. Commun.* **1**, 54-60 (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, 06/26/2023

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