

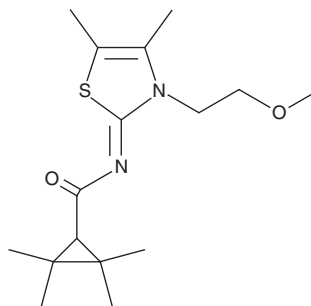
PRODUCT INFORMATION



A-836339

Item No. 13898

CAS Registry No.: 959746-77-1
Formal Name: [N(Z)]-N-[3-(2-methoxyethyl)-4,5-dimethyl-2(3H)-thiazolylidene]-2,2,3,3-tetramethyl-cyclopropanecarboxamide
MF: C₁₆H₂₆N₂O₂S
FW: 310.5
Purity: ≥98%
UV/Vis.: λ_{max}: 205, 317 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

A-836339 (Item No. 13898) is a synthetic cannabinoid (CB) which has a higher affinity for the peripheral CB₂ receptor (K_i = 0.64 nM) over the central CB₁ receptor (K_i = 270 nM).¹ It displays analgesic, anti-inflammatory, and anti-hyperalgesic effects in mice.^{1,2} At higher doses, A-836339 also decreases spontaneous locomotor activities, consistent with CB₁ activation.¹ This product is intended for forensic and research applications.

References

1. Yao, B.B., Hsieh, G., Daza, A.V., *et al.* Characterization of a cannabinoid CB₂ receptor-selective agonist, A-836339 [2,2,3,3-tetramethyl-cyclopropanecarboxylic acid [3-(2-methoxy-ethyl)-4,5-dimethyl-3H-thiazol-(2Z)-ylidene]-amide], using in vitro pharmacological assays, in vivo pain models, and pharmacological magnetic resonance imaging. *J. Pharmacol. Exp. Ther.* **328**(1), 141-151 (2009).
2. McGaraughty, S., Chu, K.L., Dart, M.J., *et al.* A CB₂ receptor agonist, A-836339, modulates wide dynamic range neuronal activity in neuropathic rats: Contributions of spinal and peripheral CB₂ receptors. *Neuroscience* **158**(4), 1652-1661 (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, 07/25/2023

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