

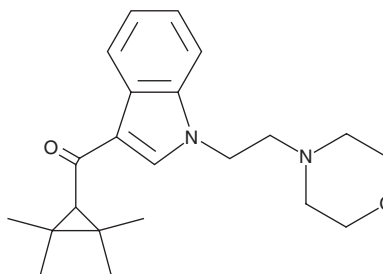
PRODUCT INFORMATION



A-796260

Item No. 11606

CAS Registry No.: 895155-26-7
Formal Name: [1-[2-(4-morpholinyl)ethyl]-1H-indol-3-yl](2,2,3,3-tetramethylcyclopropyl)methanone
Synonym: LTI-258
MF: C₂₂H₃₀N₂O₂
FW: 354.5
Purity: ≥98%
UV/Vis.: λ_{max}: 218, 246, 303 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-796260 is a synthetic cannabinoid (CB) that shows high selectivity for the CB₂ receptor (K_i = 0.77 nM) over the CB₁ receptor (K_i = 2,100 nM).¹ It also shows efficacy *in vivo* in models of inflammatory, post-operative, neuropathic, and osteoarthritic pain, activities which are selectively blocked by CB₂ receptor-selective antagonists.² This product is intended for research and forensic uses.

References

1. Poso, A. and Huffman, J.W. Targeting the cannabinoid CB₂ receptor: Modelling and structural determinants of CB₂ selective ligands. *Br. J. Pharmacol.* **153**(2), 335-346 (2008).
2. Yao, B.B., Hsieh, G.C., Frost, J.M., *et al.* *In vitro* and *in vivo* characterization of A-796260: A selective cannabinoid CB₂ receptor agonist exhibiting analgesic activity in rodent pain models. *Br. J. Pharmacol.* **153**(2), 390-401 (2008).

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, 07/18/2023

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