

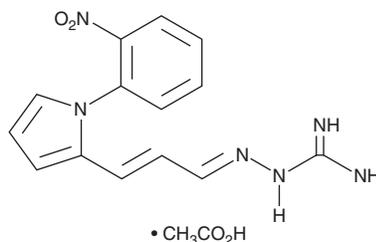
PRODUCT INFORMATION



AP1189 (acetate)

Item No. 37400

CAS Registry No.: 959850-74-9
Formal Name: 2-[3-[1-(2-nitrophenyl)-1H-pyrrol-2-yl]-2-propen-1-ylidene]-hydrazinecarboximidamide, acetate
MF: C₁₄H₁₄N₆O₂ • C₂H₄O₂
FW: 358.4
Purity: ≥95%
UV/Vis.: λ_{max}: 350 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

AP1189 (acetate) is supplied as a solid. A stock solution may be made by dissolving the AP1189 (acetate) in the solvent of choice, which should be purged with an inert gas. AP1189 (acetate) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of AP1189 (acetate) in ethanol is approximately 0.2 mg/ml and approximately 10 mg/ml in DMSO and DMF.

Description

AP1189 is a biased agonist of melanocortin receptor 1 (MC1R) and MC3R.¹ It induces ERK1 and ERK2 phosphorylation but not canonical cAMP production in HEK293 cells expressing human MC1R or MC3R when used at concentrations ranging from 10 pM to 1 mM. AP1189 (0.1-10 nM) reduces zymosan-induced IL-1β, IL-6, and TNF-α production in primary peritoneal macrophages isolated from wild-type but not *Mc3r*^{-/-} mice. It reduces zymosan-induced TNF-α production but has no effect on IL-1β or IL-6 production in primary macrophages isolated from *Mc1r*^{-/-} mice. AP1189 promotes efferocytosis in primary mouse macrophages. *In vivo*, AP1189 (1, 3, and 10 mg/kg) reduces peritoneal neutrophil infiltration in a mouse model of zymosan-induced acute peritonitis. AP1189 (50 mg/kg) reduces paw volume, the number of paws affected, and disease severity in a mouse model of arthritogenic serum-induced arthritis.

Reference

1. Montero-Melendez, T., Gobbetti, T., Cooray, S.N., *et al.* Biased agonism as a novel strategy to harness the proresolving properties of melanocortin receptors without eliciting melanogenic effects. *J. Immunol.* **194**(7), 3381-3388 (2015).

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, 06/15/2023

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