

PRODUCT INFORMATION

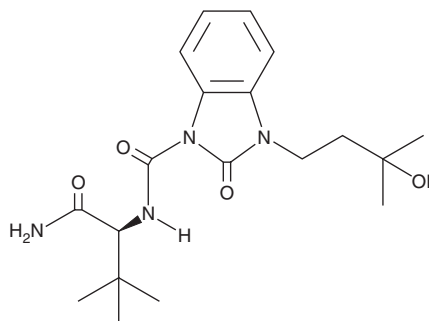


PF-03550096

Item No. 15620

CAS Registry No.: 910376-39-5
Formal Name: N-[(1S)-1-(aminocarbonyl)-2,2-dimethylpropyl]-2,3-dihydro-3-(3-hydroxy-3-methylbutyl)-2-oxo-1H-benzimidazole-1-carboxamide

MF: C₁₉H₂₈N₄O₄
FW: 376.5
Purity: ≥98%
UV/Vis.: λ_{max}: 212 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

PF-03550096 is a synthetic cannabinoid (CB) that selectively targets the peripheral CB₂ receptor over the central CB₁ receptor (K_is = 7.9 and 1,500 nM for human isoforms, respectively).^{1,2} It shows similar selectivity for inhibiting forskolin-stimulated cAMP production in CHO cells expressing human CB₁ and CB₂ receptors (EC₅₀s = 450 and 2.1 nM, respectively).² PF-03550096 is orally bioavailable and significantly reduces colonic pain induced by 2,4,6-trinitrobenzene sulfonic acid in rats.² This product is intended for forensic and research applications.

References

1. Ando, K., Buchler, I., Hedge, S.G., *et al.* Benzimidazolone derivatives. *Pfizer Products Inc.* **WO/2008/032164** (2009).
2. Kikuchi, A., Ohashi, K., Sugie, Y., *et al.* Pharmacological evaluation of a novel cannabinoid 2 (CB₂) ligand, PF-03550096, in vitro and in vivo by using a rat model of visceral hypersensitivity. *J. Pharmacol. Sci.* **106(2)**, 219-224 (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, 06/28/2023

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