

PRODUCT INFORMATION

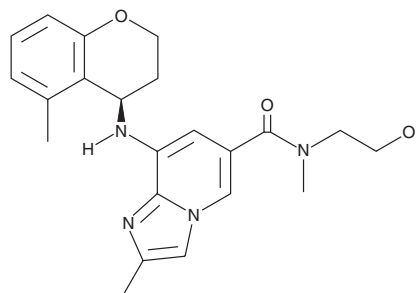


PF-3716556

Item No. 17710

CAS Registry No.: 928774-43-0
Formal Name: 8-[[[(4R)-3,4-dihydro-5-methyl-2H-1-benzopyran-4-yl]amino]-N-(2-hydroxyethyl)-N,2-dimethylimidazo[1,2-a]pyridine-6-carboxamide

MF: C₂₂H₂₆N₄O₃
FW: 394.5
Purity: ≥98%
UV/Vis.: λ_{max}: 254, 297 nm
Supplied as: A crystalline 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

PF-3716556 is supplied as a crystalline solid. A stock solution may be made by dissolving the PF-3716556 in the solvent of choice, which should be purged with an inert gas. PF-3716556 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of PF-3716556 in ethanol is approximately 10 mg/ml and approximately 30 mg/ml in DMSO and DMF.

PF-3716556 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, PF-3716556 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. PF-3716556 has a solubility of approximately 0.04 mg/ml in a 1:20 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

PF-3716556 is a potent and selective hydrogen-potassium pump (H⁺/K⁺-ATPase) inhibitor in a new class of acid suppressants, the potassium-competitive acid blockers or acid pump antagonists.¹ PF-3716556 inhibited H⁺/K⁺-ATPase activity in porcine membrane vesicles, as well as canine and human recombinant cells under ion-leaky conditions (pI_{C₅₀ = 6.026, 6.038, and 6.009, respectively). PF-03716556 showed 3-fold greater inhibition than revaprazan in ion-tight (neutral pH) conditions.}

Reference

1. Mori, H., Tonai-Kachi, H., Ochi, Y., *et al.* N-(2-hydroxyethyl)-N,2-dimethyl-8-[[[(4R)-5-methyl-3,4-dihydro-2H-chromen-4-yl]amino]imidazo[1,2-a]pyridine-6-carboxamide (PF-03716556), a novel, potent, and selective acid pump antagonist for the treatment of gastroesophageal reflux disease. *J. Pharmacol. Exp. Ther.* **328**(2), 671-679 (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, 10/06/2022

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