

PRODUCT INFORMATION



A-1596586

Item No. 41316

CAS Registry No.: 1918142-34-3
Formal Name: 1-(2,2-difluoro-1,3-benzodioxol-5-yl)-N-[(2R,4R)-2-(3,4-dimethoxyphenyl)-3,4-dihydro-7-methoxy-2H-1-benzopyran-4-yl]-cyclopropanecarboxamide

MF: C₂₉H₂₇F₂NO₇

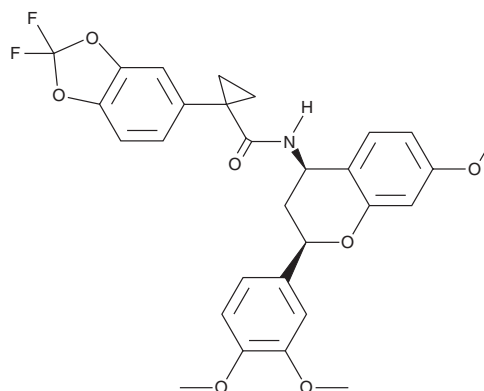
FW: 539.5

Purity: ≥98%

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

A-1596586 is supplied as a solid. A stock solution may be made by dissolving the A-1596586 in the solvent of choice, which should be purged with an inert gas. A-1596586 is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

A-1596586 is a corrector of the cystic fibrosis transmembrane conductance regulator (CFTR) bearing the F508 deletion (F508del) mutation.¹ It increases CFTR levels in CFBE41o- bronchial epithelial cells expressing F508del-CFTR fused to HRP (EC₅₀ = 0.13 μM) and the transepithelial current in primary human bronchial epithelial cells isolated from a patient with cystic fibrosis with the F508del-CFTR mutation (EC₅₀ = 0.028 μM).

Reference

1. Wang, X., Liu, B., Searle, X., et al. Discovery of 4-[(2R,4R)-4-({[1-(2,2-difluoro-1,3-benzodioxol-5-yl)cyclopropyl]carbonyl}amino)-7-(difluoromethoxy)-3,4-dihydro-2H-chromen-2-yl]benzoic acid (ABBV/GLPG-2222), a potent cystic fibrosis transmembrane conductance regulator (CFTR) corrector for the treatment of cystic fibrosis. *J. Med. Chem.* **61**(4), 1436-1449 (2017).

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, 08/30/2024

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