

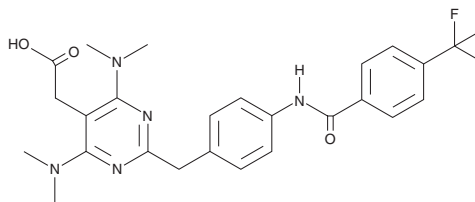
PRODUCT INFORMATION



BI-671800

Item No. 27638

CAS Registry No.: 1093108-50-9
Formal Name: 4,6-bis(dimethylamino)-2-[[4-[[4-(trifluoromethyl)benzoyl]amino]phenyl]methyl]-5-pyrimidineacetic acid
Synonym: AP-761
MF: C₂₅H₂₆F₃N₅O₃
FW: 501.5
Purity: ≥98%
UV/Vis.: λ_{max}: 251, 279 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

BI-671800 is supplied as a crystalline solid. A stock solution may be made by dissolving the BI-671800 in the solvent of choice, which should be purged with an inert gas. BI-671800 is soluble in the organic solvent DMSO.

Description

BI-671800 is an antagonist of the prostaglandin D₂ (PGD₂; Item No. 12010) receptor CRTH₂/DP₂ (IC₅₀ = 42 nM in a radioligand binding assay).¹ It binds selectively to CRTH₂/DP₂ over DP₁, thromboxane, and prostacyclin receptors (IC₅₀s = 25.6, 62.6, and >100 μM, respectively). BI-671800 inhibits PGD₂-induced shape change in human eosinophils (IC₅₀ = 744 nM).

Reference

1. Stearns, B.A., Baccei, C., Bain, G., *et al.* Novel tricyclic antagonists of the prostaglandin D₂ receptor DP₂ with efficacy in a murine model of allergic rhinitis. *Bioorg. Med. Chem. Lett.* **19(16)**, 4647-4651 (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, 12/22/2022

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