

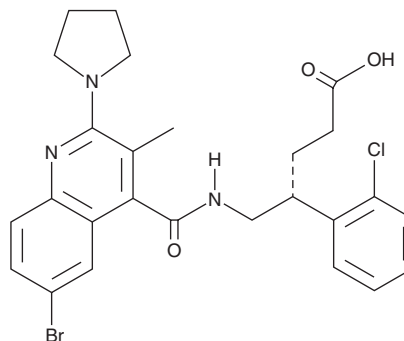
PRODUCT INFORMATION



BAY-6672

Item No. 42731

CAS Registry No.: 2247517-53-7
Formal Name: (+)-γ-[[[6-bromo-3-methyl-2-(1-pyrrolidinyl)-4-quinolyl]carbonyl]amino]methyl]-2-chloro-benzenebutanoic acid
MF: C₂₆H₂₇BrClN₃O₃
FW: 544.9
Purity: ≥95%
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

BAY-6672 is supplied as a solid. A stock solution may be made by dissolving the BAY-6672 in the solvent of choice, which should be purged with an inert gas. BAY-6672 is slightly soluble (0.1-1 mg/ml) in acetonitrile and DMSO.

Description

BAY-6672 is an FP receptor antagonist (IC₅₀ = 11 nM).¹ It is selective for the FP receptor over the prostaglandin E₂ (PGE₂) receptor subtypes EP₁, EP₂, EP₃, and EP₄ (IC₅₀s = >10, >10, >10, and >9.4 μM, respectively), as well as IP and DP receptors (IC₅₀s = >10 μM for both). *In vivo*, BAY-6672 (30 mg/kg) reduces lung IL-1β levels and fibrosis in a mouse model of silica-induced pulmonary fibrosis.

Reference

1. Beck, H., Thaler, T., Meibom, D., *et al.* Potent and selective human prostaglandin F (FP) receptor antagonist (BAY-6672) for the treatment of idiopathic pulmonary fibrosis (IPF). *J. Med. Chem.* **63**(20), 11639-11662 (2020).

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, 04/02/2025

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