

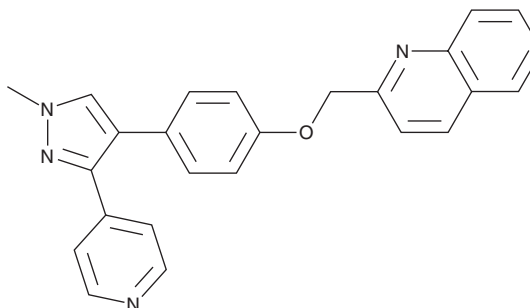
PRODUCT INFORMATION



PF-02545920

Item No. 42047

CAS Registry No.: 898563-49-0
Formal Name: 2-[[4-[1-methyl-3-(4-pyridinyl)-1H-pyrazol-4-yl]phenoxy]methyl]-quinoline
MF: C₂₅H₂₀N₄O
FW: 392.5
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

PF-02545920 is supplied as a solid. A stock solution may be made by dissolving the PF-02545920 in the solvent of choice, which should be purged with an inert gas. PF-02545920 is sparingly soluble (1-10 mg/ml) in DMSO.

Description

PF-02545920 is an inhibitor of phosphodiesterase 10A (PDE10A; IC₅₀ = 29 nM).¹ It inhibits PCP-induced hyperactivity in mice when administered at doses of 30 or 100 mg/kg. PF-02545920 (0.1 or 0.3 mg/kg) reverses PCP-induced deficits in the novel object recognition test in neonatal mice.

Reference

1. Hamaguchi, W., Masuda, N., Samizu, K., *et al.* Synthesis and *in vivo* evaluation of novel quinoline derivatives as phosphodiesterase 10A inhibitors. *Chem. Pharm. Bull. (Tokyo)* **62(12)**, 1200-1213 (2014).

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, 01/21/2025

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