

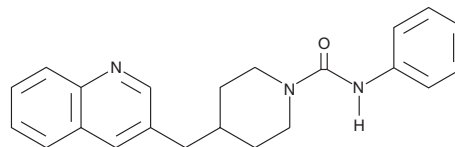
PRODUCT INFORMATION



PF-750

Item No. 10010908

CAS Registry No.: 959151-50-9
Formal Name: N-phenyl-4-(quinolin-2-ylmethyl)piperidine-1-carboxamide
MF: C₂₂H₂₃N₃O
FW: 345.4
Purity: ≥98%
UV/Vis.: λ_{max}: 206, 234 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-750 is supplied as a crystalline solid. A stock solution may be made by dissolving the PF-750 in the solvent of choice, which should be purged with an inert gas. PF-750 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of PF-750 in ethanol is approximately 1 mg/ml and approximately 20 mg/ml in DMSO and DMF.

PF-750 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, PF-750 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. PF-750 has a solubility of approximately 0.125 mg/ml in a 1:5 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Fatty acid amide hydrolase (FAAH) is the enzyme responsible for hydrolysis and inactivation of fatty acid amides including anandamide and oleamide. PF-750 is a potent, time-dependent, irreversible FAAH inhibitor with IC₅₀ values of 0.6 and 0.016 μM when preincubated with recombinant human FAAH for 5 and 60 minutes, respectively.¹ Activity-based profiling of various human and murine tissue proteome samples revealed that PF-750 is highly selective for FAAH relative to other serine hydrolases, showing no discernable off-site activity up to 500 μM.¹

Reference

1. Ahn, K., Johnson, D.S., Fitzgerald, L.R., *et al.* Novel mechanistic class of fatty acid amide hydrolase inhibitors with remarkable selectivity. *Biochemistry* **46**, 13019-13030 (2007).

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, 05/15/2024

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