

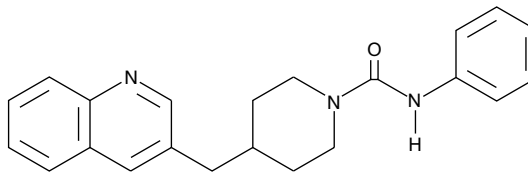
Product Information



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%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid
UV/Vis.: λ_{max}: 206, 234 nm



Laboratory Procedures

For long term storage, we suggest that PF-750 be stored as supplied at -20°C. It should be stable for at least two years. PF-750 is supplied as a crystalline solid. A stock solution may be made by dissolving the PF-750 in an organic solvent 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.

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).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/10010908

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will meet our specifications at the time of delivery.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a refund of the purchase price, or at Cayman's option, the replacement, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy** located on our website and in our catalog.

Copyright Cayman Chemical Company, 02/16/2012