

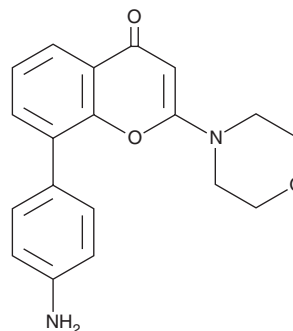
PRODUCT INFORMATION



PI-828

Item No. 21811

CAS Registry No.: 942289-87-4
Formal Name: 8-(4-aminophenyl)-2-(4-morpholinyl)-4H-1-benzopyran-4-one
MF: C₁₉H₁₈N₂O₃
FW: 322.4
Purity: ≥98%
UV/Vis.: λ_{max}: 245, 305 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

PI-828 is supplied as a crystalline solid. A stock solution may be made by dissolving the PI-828 in the solvent of choice, which should be purged with an inert gas. PI-828 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of PI-828 in these solvents is approximately 25 mg/ml. PI-828 is also slightly soluble in ethanol.

Description

PI-828 is a potent inhibitor of phosphatidylinositol 3-kinase (PI3K; IC₅₀s = 9.8, 183, 227, and 1,967 nM for p110β, p110α, p110δ, and p110γ, respectively).¹ This molecule has been immobilized on epoxy-containing solid surface via the 4'amine for binding studies and shown to interact with non-PI3K cellular proteins such as mTOR, ALDH1, BRD4, CKIIα, and GSK-3β.

Reference

1. Gharbi, S.I., Zvelebil, M.J., Shuttleworth, S.J., *et al.* Exploring the specificity of the PI3K family inhibitor LY294002. *Biochem. J.* **404**(1), 15-21 (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, 12/05/2022

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