

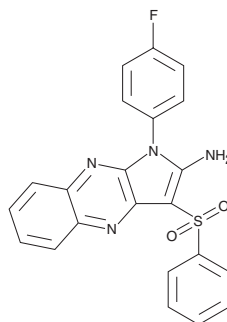
Product Information



CAY10602

Catalog No. 10009796

CAS Registry No.: 374922-43-7
Formal Name: 1-(4-fluorophenyl)-3-(phenylsulfonyl)-1H-pyrrolo[2,3-b]quinoxalin-2-amine
MF: C₂₂H₁₅FN₄O₂S
FW: 418.4
Purity: ≥95%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid



Laboratory Procedures

For long term storage, we suggest that CAY10602 be stored as supplied at -20°C. It should be stable for at least two years.

CAY10602 is supplied as a crystalline solid. A stock solution may be made by dissolving the CAY10602 in an organic solvent purged with an inert gas. CAY10602 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of CAY10602 in these solvents is approximately 20 mg/ml.

CAY10602 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CAY10602 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. CAY10602 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Sirtuin 1 (SIRT1) is an evolutionarily conserved NAD-dependent protein deacetylase that has been implicated in cell metabolism, differentiation, stress and DNA damage response, and the control of multidrug resistance in cancer. SIRT1 alters gene expression by deacetylating nuclear histones, NF-κB, p53, and FOXO transcription factors. CAY10602 was derived from high throughput screening for compounds that increase SIRT1-mediated deacetylation of a SIRT1-specific substrate.¹ Functional assays show that CAY10602 dose-dependently suppresses the NF-κB-dependent induction of TNF-α by lipopolysaccharide in THP-1 cells, with approximately 75% inhibition achieved at 60 μM, without cytotoxicity.¹ CAY10602 is at least 10 times more potent as a suppressor of TNF-α release than resveratrol. The effects of CAY10602 in other cellular processes remain to be determined.

Reference

1. Nayagam, V.M., Wang, X., Tan, Y.C., *et al.* SIRT1 modulating compounds from high-throughput screening as anti-inflammatory and insulin-sensitizing agents. *Journal of Biomolecular Screening* **11**(8), 959-967 (2006).

Related Products

trans-trismethoxy Resveratrol - Cat. No. 10188 • *trans*-trismethoxy Resveratrol-d₄ - Cat. No. 13129 • *cis*-trismethoxy Resveratrol - Cat. No. 13199 • *trans*-Resveratrol - Cat. No. 70675 • HDAC Activity/Inhibitor Screening Assay Kit - Cat. No. 789701 • *cis*-Resveratrol - Cat. No. 10004235 • CAY10591 - Cat. No. 10009797 • SIRT1 Direct Fluorescent Screening Assay Kit - Cat. No. 10010401 • SIRT1 FRET-Based Screening Assay Kit - Cat. No. 10010991 • SIRT1 (human recombinant) - Cat. No. 10011190

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, 10/13/2009

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