

Product Information



Caylin-2

Item No. 10005002

Formal Name: 4-[4,5-bis(4-trifluoromethyl-phenyl)-2-(2-isopropoxy-4-methoxy-phenyl)-4,5-dihydro-imidazole-1-carboxyl]-piperazin-2-one

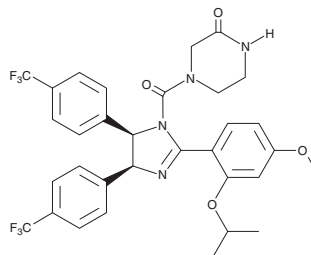
MF: C₃₂H₃₀F₆N₄O₄

FW: 648.6

Purity: ≥98%

Stability: ≥1 year at -20°C

Supplied as: A crystalline solid



Laboratory Procedures

For long term storage, we suggest that caylin-2 be stored as supplied at -20°C. It will be stable for at least one year.

Caylin-2 is supplied as a crystalline solid. A stock solution may be made by dissolving the caylin-2 in an organic solvent purged with an inert gas. Caylin-2 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of caylin-2 in these solvents is at least 30 mg/ml.

Caylin-2 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, caylin-2 should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Caylin-2 has a solubility of 0.025 mg/ml in a 1:3 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Nutlin-3 is an activator of p53 that functions by inhibiting the interaction of p53 with MDM2, a negative regulator of p53 activity.¹ Caylin-2 is a nutlin-3 analog in which trifluoromethyl groups have been substituted for chlorine on two of the phenyl rings. At high concentrations, caylin-2 inhibits the growth of HCT-116 cells with an IC₅₀ of approximately 8 μM, making it about 10-fold less potent than nutlin-3. Interestingly, at concentrations between 5-100 nM, caylin-2 promotes the growth of HCT-116 cells approximately 40% compared to untreated cells.² The mechanism of the growth promoting properties of caylin-2 have not yet been elucidated.

References

1. Vassilev, L.T., Vu, B.T., Graves, B., *et al.* *In vivo* activation of the p53 pathway by small-molecule antagonists of MDM2. *Science* **303**, 844-848 (2004).
2. Meade, E.A., [Unpublished] (2004).

Related Products

p53 (human recombinant) - Item No. 10357 • PRIMA-1 - Item No. 63520 • Hepsin Polyclonal Antibody - Affinity-Purified (aa 241-260) - Item No. 100022 • β-Catenin Polyclonal Antibody - Item No. 100029

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 catalog.

Copyright Cayman Chemical Company, 03/03/2011