

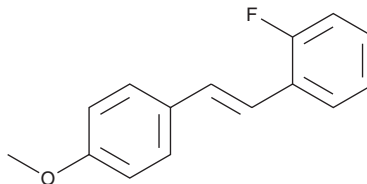
Product Information



CAY10512

Item No. 10009536

CAS Registry No.: 139141-12-1
Formal Name: 1-fluoro-2-[2-(4-methoxyphenyl)ethenyl]-benzene
MF: C₁₅H₁₃FO
FW: 228.3
Purity: ≥97%
Stability: ≥2 years at -20°C
Supplied as: A crystalline solid



Laboratory Procedures

For long term storage, we suggest that CAY10512 be stored as supplied at -20°C. It will be stable for at least two years.

CAY10512 is supplied as a crystalline solid. A stock solution may be made by dissolving the CAY10512 in an organic solvent purged with an inert gas. CAY10512 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of CAY10512 in ethanol, DMSO, and DMF is approximately 0.5, 2, and 3 mg/ml, respectively.

CAY10512 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CAY10512 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. CAY10512 has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

The nuclear factor-κB (NF-κB) regulates the expression of numerous genes involved in immunity and inflammation, cellular stress responses, growth, and apoptosis. Resveratrol, a polyphenolic *trans*-stilbene, is a known inhibitor of the activation of NF-κB and exhibits activity against a wide variety of cancer cells. CAY10512 is a substituted *trans*-stilbene analog of resveratrol that is 100-fold more potent as measured by antioxidant activity. The IC₅₀ value for inhibition of TNF-α-induced activation of NF-κB by CAY10512 is 0.15 μM compared to 20 μM by resveratrol.¹

Reference

1. Heynekamp, J.J., Weber, W.M., Hunsaker, L.A., *et al.* Substituted *trans*-stilbenes, including analogues of the natural product resveratrol, inhibit the human tumor necrosis factor α-induced activation of transcription factor nuclear factor κB. *J. Med. Chem.* **49**, 7182-7189 (2006).

Related Product

Trismethoxy Resveratrol - Item No. 10188 • Resveratrol - Catalog No. 70675 • PDM 11 - Item No. 10006341 • PDM 2 - Item No. 10006342 • CAY10465 - Item No. 10006546

WARNING: THIS PRODUCT IS NOT FOR HUMAN OR ANIMAL DISEASE DIAGNOSIS OR THERAPEUTIC DRUG 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 under separate cover to the MSDS supervisor at 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, 09/30/2011

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