

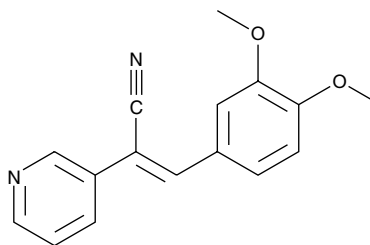
Product Information



RG-13022

Item No. 10010309

CAS Registry No.: 149286-90-8
Formal Name: α Z-[(3,4-dimethoxyphenyl)methylene]-3-pyridineacetonitrile
MF: C₁₆H₁₄N₂O₂
FW: 266.3
Purity: \geq 98%
Stability: \geq 2 years at -20°C
Supplied as: A crystalline solid
UV/Vis.: λ_{max} : 248, 356 nm



Laboratory Procedures

For long term storage, we suggest that RG-13022 be stored as supplied at -20°C. It will be stable for at least two years.

RG-13022 is supplied as a crystalline solid. A stock solution may be made by dissolving the RG-13022 in an organic solvent purged with an inert gas. RG-13022 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of RG-13022 in ethanol is approximately 2 mg/ml and approximately 10 mg/ml in DMSO and DMF.

RG-13022 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, RG-13022 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. RG-13022 has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Protein tyrosine kinase (PTK) inhibitors are potential antiproliferative agents for diseases caused by the hyperactivity of PTKs. Tyrphostins are a class of antiproliferative compounds that act as PTK blockers. PTK inhibitors that preferentially inhibit the epidermal growth factor (EGF) receptor kinase block EGF-dependent cell proliferation.¹ RG-13022 is an inhibitor of EGF receptor kinase with an IC₅₀ value of 1 μ M in HT-22 cells.²

References

1. Gazit, A., Yaish, P., Gilon, C., *et al.* Tyrphostins I: Synthesis and biological activity of protein tyrosine kinase inhibitors. *J. Med. Chem.* **32(10)**, 2344-2352 (1989).
2. Sagara, Y., Ishige, K., Tsai, C., *et al.* Tyrphostins protect neuronal cells from oxidative stress. *J. Biol. Chem.* **277(39)**, 36204-36215 (2002).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/10010309

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, 04/03/2012

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