

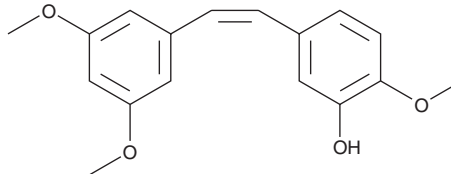
PRODUCT INFORMATION



CAY10616

Item No. 13291

CAS Registry No.: 586410-08-4
Formal Name: 5-[(1Z)-2-(3,5-dimethoxyphenyl)ethenyl]-2-methoxy-phenol
MF: C₁₇H₁₈O₄
FW: 286.3
Purity: ≥98%
UV/Vis.: λ_{max}: 291 nm
Supplied as: A solution in methyl acetate
Storage: -20°C
Stability: ≥1 year



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

CAY10616 is supplied as a solution in methyl acetate. To change the solvent, simply evaporate the methyl acetate under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide (DMF) purged with an inert gas can be used. The solubility of CAY10616 in these solvents is approximately 10, 14, and 20 mg/ml, respectively.

CAY10616 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the methyl acetate solution of CAY10616 should be diluted with the aqueous buffer of choice. CAY10616 has a solubility of 0.1 mg/ml in a 1:9 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Resveratrol is a natural polyphenolic antioxidant that has anti-cancer properties. CAY10616 is an analog of resveratrol which potently induces apoptosis in promyelocytic leukemia HL-60 cells (IC₅₀ = 40 nM, vs. 50 μM for resveratrol).¹ In addition, CAY10599 induces apoptosis (IC₅₀ = 30 nM) in HL-60R, a cell line derived from HL-60 that expresses a multidrug-resistance phenotype (resistant to daunorubicin, etoposide, and citarabine).¹

Reference

1. Roberti, M., Pizzirani, D., Simoni, D., *et al.* Synthesis and biological evaluation of resveratrol and analogues as apoptosis-inducing agents. *J. Med. Chem.* **46**(16), 3546-3554 (2003).

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, 04/26/2023

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