

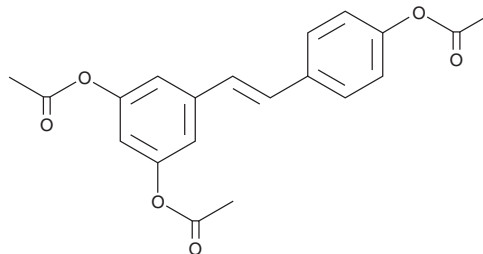
# PRODUCT INFORMATION



## Acetyl Resveratrol

Item No. 26379

**CAS Registry No.:** 42206-94-0  
**Formal Name:** 5-[(1E)-2-[4-(acetyloxy)phenyl]ethenyl]-1,3-benzenediol, 1,3-diacetate  
**Synonyms:** Resveratrol triacetate, Triacetylresveratrol, 3,5,4'-Tri-O-acetylresveratrol  
**MF:** C<sub>20</sub>H<sub>18</sub>O<sub>6</sub>  
**FW:** 354.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 224, 300, 310 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years



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

### Laboratory Procedures

Acetyl resveratrol is supplied as a crystalline solid. A stock solution may be made by dissolving the acetyl resveratrol in the solvent of choice. Acetyl resveratrol is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of acetyl resveratrol in these solvents is approximately 50 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of acetyl resveratrol can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of acetyl resveratrol in PBS, pH 7.2, is >0.1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Acetyl resveratrol is a triacetylated derivative of resveratrol that has diverse biological activities, including anticancer and antioxidant properties.<sup>1-3</sup> It inhibits the growth of LNCaP, DU145, and PC3M prostate cancer cells (IC<sub>50</sub>s = 10.5, 14.2, and 26.8 μM, respectively).<sup>2</sup> Acetyl resveratrol induces cytotoxicity in ALL-5 acute lymphoblastic leukemia cells (IC<sub>50</sub> = 3.4 μM) and induces cell cycle arrest at the G<sub>1</sub> phase when used at a concentration of 17 μM.<sup>1</sup> It also increases survival in mice exposed to total body γ-irradiation when administered at a dose of 10 mg/kg prior to irradiation.<sup>3</sup>

### References

1. Urbaniak, A., Delgado, M., Kacprzak, K., *et al.* Activity of resveratrol triesters against primary acute lymphoblastic leukemia cells. *Bioor. Med. Chem. Lett.* **27(12)**, 2766-2770 (2017).
2. Dias, S.J., Li, K., Rimando, A.M., *et al.* Trimethoxy-resveratrol and piceatannol administered orally suppress and inhibit tumor formation and growth in prostate cancer xenografts. *Prostate* **73(11)**, 1135-1146 (2016).
3. Koide, K., Osman, S., Garner, A.L., *et al.* The use of 3,5,4'-Tri-O-acetylresveratrol as a potential pro-drug for resveratrol protects mice from γ-irradiation-induced death. *ACS Med. Chem. Lett.* **2(4)**, 270-274 (2011).

#### 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, 12/19/2022

#### 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