

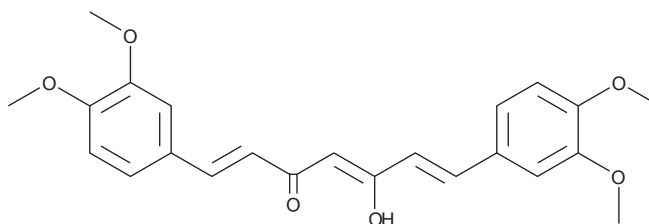
# Product Information



## ASC-J9

Catalog No. 10009986

**CAS Registry No.:** 917813-54-8  
**Formal Name:** 1,7-bis(3,4-dimethoxyphenyl)-5-hydroxy-1E,4E,6E-heptatrien-3-one  
**MF:** C<sub>23</sub>H<sub>24</sub>O<sub>6</sub>  
**FW:** 396.4  
**Purity:** ≥95%  
**Stability:** ≥2 years at -20°C  
**Supplied as:** A crystalline solid



### Laboratory Procedures

For long term storage, we suggest that ASC-J9 be stored as supplied at -20°C. It will be stable for at least two years.

ASC-J9 is supplied as a crystalline solid. A stock solution may be made by dissolving the ASC-J9 in an organic solvent purged with an inert gas. ASC-J9 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of ASC-J9 in these solvents is approximately 0.25, 5, and 10 mg/ml respectively.

ASC-J9 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, ASC-J9 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. ASC-J9 has a solubility of approximately 0.30 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.

Naturally occurring phytochemicals such as turmeric (curcumin) have been found to inhibit the growth of tumor cells. ASC-J9 is an analog of curcumin obtained by methylation of both free phenolic groups in the parent compound. A 30-fold increase in potency was observed in the growth suppression of HCT116 tumor cells following this modification.<sup>1</sup>

### Reference

1. Otori, H., Yamakoshi, H., Tomizawa, M., *et al.* Synthesis and biological analysis of new curcumin analogues bearing an enhanced potential for the medicinal treatment of cancer. *Mol. Cancer Therapeutics* **5(10)**, 2563-2571 (2006).

### Related Products

Celastrol - Cat. No. 70950 • Curcumin - Cat. No. 81025 • Curcumin (technical grade) - Cat. No. 81025.1

## 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 NOT INTENDED OR APPROVED FOR HUMAN OR VETERINARY USE. USE OF THIS PRODUCT FOR HUMAN OR ANIMAL TESTING IS EXTREMELY HAZARDOUS AND MAY RESULT IN DISEASE, SEVERE INJURY, OR DEATH.**

#### 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, 08/14/2007