

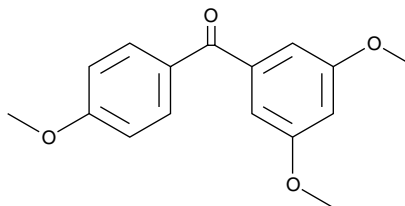
Product Information



3,4',5-Trimethoxybenzophenone

Item No. 10004185

CAS Registry No.: 94709-12-3
Formal Name: (3,5-dimethoxyphenyl)(4-methoxyphenyl)-methanone
MF: C₁₆H₁₆O₄
FW: 272.3
Purity: ≥95%
Stability: ≥1 year at -20°C
Supplied as: A crystalline solid
UV/Vis.: λ_{max}: 222, 384 nm



Laboratory Procedures

For long term storage, we suggest that 3,4',5-trimethoxybenzophenone be stored as supplied at -20°C. It should be stable for at least one year.

3,4',5-Trimethoxybenzophenone is supplied as a crystalline solid. A stock solution may be made by dissolving the 3,4',5-trimethoxybenzophenone in an organic solvent purged with an inert gas. 3,4',5-Trimethoxybenzophenone is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of 3,4',5-trimethoxybenzophenone in these solvents is at least 20 mg/ml.

3,4',5-Trimethoxybenzophenone is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, 3,4',5-trimethoxybenzophenone should first be dissolved in DMF and then diluted with the aqueous buffer of choice. 3,4',5-Trimethoxybenzophenone has a solubility of 0.1 mg/ml in a 1:3 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Resveratrol is a potent phenolic antioxidant found in natural sources that has antiproliferative activity.¹ When the three phenolic hydroxyl groups of resveratrol are converted to methyl ethers, the inhibition of cell growth is enhanced.² 3,4',5-Trimethoxybenzophenone is an analog of trimethoxy resveratrol. It inhibits the growth of a variety of human tumor cell lines at concentrations from 0.4 to 2 μg/ml, which is 5-6 times more potent than resveratrol.³

References

1. Rotondo, S., Rajtar, G., Manarini, S., *et al.* Effect of *trans*-resveratrol, a natural polyphenolic compound, on human polymorphonuclear leukocyte function. *Br. J. Pharmacol.* **123**, 1691-1699 (1998).
2. Nam, K.A., Kim, S., Heo, Y.H., *et al.* Resveratrol analog, 3,5,2',4'-tetramethoxy-*trans*-stilbene, potentiates the inhibition of cell growth and induces apoptosis in human cancer cells. *Arch. Pharm. Res.* **24(5)**, 441-445 (2001).
3. Pettit, G.R., Grealish, M.P., Jung, M.K., *et al.* Antineoplastic agents. 465. Structural modification of resveratrol: Sodium resverastatin phosphate. *J. Med. Chem.* **45**, 2534-2542 (2002).

Related Products

TMS - Item No. 10038 • Trimethoxy Resveratrol - Item No. 10188 • Resveratrol - Item No. 70675

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 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, 06/20/2011