

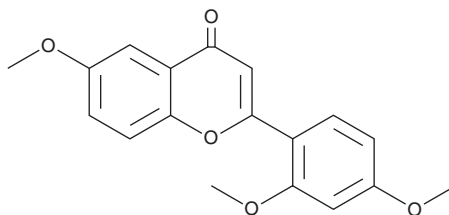
PRODUCT INFORMATION



6,2',4'-Trimethoxyflavone

Item No. 33150

CAS Registry No.: 720675-74-1
Formal Name: 2-(2,4-dimethoxyphenyl)-6-methoxy-4H-1-benzopyran-4-one
MF: C₁₈H₁₆O₅
FW: 312.3
Purity: ≥95%
UV/Vis.: λ_{max}: 226, 271, 334 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Synthetic



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

Laboratory Procedures

6,2',4'-Trimethoxyflavone is supplied as a crystalline solid. A stock solution may be made by dissolving the 6,2',4'-trimethoxyflavone in the solvent of choice, which should be purged with an inert gas. 6,2',4'-Trimethoxyflavone is soluble in organic solvents chloroform, ethanol, DMSO and dimethyl formamide (DMF). The solubility of 6,2',4'-trimethoxyflavone in chloroform is approximately 10 mg/ml. 6,2',4'-Trimethoxyflavone is slightly soluble in ethanol, DMSO, and DMF.

Description

6,2',4'-Trimethoxyflavone is a flavonoid that has been found in *T. acutifolius* and has diverse biological activities.¹⁻³ It acts as an antagonist of the aryl hydrocarbon receptor (AhR), inhibiting benzo[a]pyrene-induced AhR-dependent transcription in a reporter assay when used at concentrations of 2 and 10 μM.¹ 6,2',4'-Trimethoxyflavone decreases LPS-induced production of TNF-α in THP-1 cells (IC₅₀ = 47.7 μg/ml).² It inhibits cell migration and invasion, but not proliferation, of HN-30 head and neck squamous cell carcinoma (HNSCC) cells *in vitro* when used at a concentration of 10 μM.³

References

1. Murray, I.A., Flaveny, C.A., DiNatale, B.C., *et al* Antagonism of aryl hydrocarbon receptor signaling by 6,2',4'-trimethoxyflavone. *J. Pharmacol. Exp. Ther.* **332**(1), 135-144 (2010).
2. Apaza T, L., Serban, A.M., Cabanillas, A.H., *et al*. Flavonoids of *Tripodanthus acutifolius* inhibit TNF-α production in LPS-activated THP-1 and B16-F10 cells. *J. Ethnopharmacol.* **242**, 112036 (2019).
3. DiNatale, B.C., Smith, K., John, K., *et al*. Ah receptor antagonism represses head and neck tumor cell aggressive phenotype. *Mol. Cancer Res.* **10**(10), 1369-1379 (2012).

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, 10/11/2022

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