

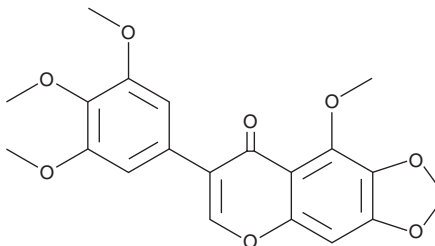
PRODUCT INFORMATION



Irisflorentin

Item No. 46194

CAS Registry No.: 41743-73-1
Formal Name: 9-methoxy-7-(3,4,5-trimethoxyphenyl)-8H-1,3-dioxolo[4,5-g][1]benzopyran-8-one
Synonyms: 41743-73-1
MF: C₂₀H₁₈O₈
FW: 386.4
Purity: ≥98%
Supplied as: A 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

Irisflorentin is supplied as a solid. A stock solution may be made by dissolving the irisflorentin in the solvent of choice, which should be purged with an inert gas. Irisflorentin is sparingly soluble (1-10 mg/ml) in DMSO.

Description

Irisflorentin is an isoflavone that has been found in *B. chinensis* and has anti-inflammatory and cytoprotective activities.¹⁻³ It reduces LPS-induced increases in the levels of nitric oxide (NO), inducible NO synthase (iNOS), TNF- α , IL-1 β , and IL-6 in RAW 264.7 macrophages at 10, 20, and 40 μ M.¹ Irisflorentin increases phagocytosis of *E. coli* by RAW 264.7 macrophages at 40 μ M and prevents M1 polarization in mouse bone marrow-derived macrophages (BMDMs).² It also reduces the secretion of IL-6, TNF- α , and IL-10 induced by heat-inactivated methicillin-resistant *S. aureus* (HK-MRSA) in mouse BMDMs. Irisflorentin activates the PI3K, mTOR, and BDNF signaling pathways, increases cell viability, and reduces lactate dehydrogenase (LDH) release, reactive oxygen species (ROS) production, and apoptosis in a PC12 cellular model of ischemic stroke using oxygen-glucose deprivation and reoxygenation.³

References

1. Gao, Y., Fang, L., Liu, F., et al. Suppressive effects of irisflorentin on LPS-induced inflammatory responses in RAW 264.7 macrophages. *Exp. Biol. Med. (Maywood)* **239**(8), 1018-1024 (2014).
2. Xiang, T., Zou, Y., Jiang, X., et al. Irisflorentin promotes bacterial phagocytosis and inhibits inflammatory responses in macrophages during bacterial infection. *Heliyon* **10**(1), e23225 (2023).
3. Y., Z., Lu, T., Luo, X., et al. Elucidating the mechanism of anti-ischemic stroke action of Irisflorentin: A combined approach using network pharmacology, molecular docking, and in vitro validation. *Eur. J. Med. Chem. Rep.* **16**, 1000318 (2026).

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, 08/05/2026

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