

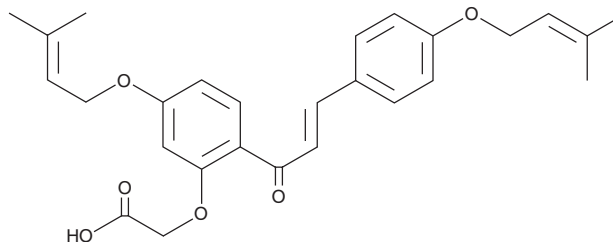
PRODUCT INFORMATION



Sofalcone

Item No. 11882

CAS Registry No.: 64506-49-6
Formal Name: 2-[5-[(3-methyl-2-buten-1-yl)oxy]-2-[3-[4-[(3-methyl-2-buten-1-yl)oxy]phenyl]-1-oxo-2-propen-1-yl]phenoxy]-acetic acid
Synonyms: Solon, SU 88
MF: C₂₇H₃₀O₆
FW: 450.5
Purity: ≥98%
UV/Vis.: λ_{max}: 238, 348 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

Sofalcone is supplied as a crystalline solid. A stock solution may be made by dissolving the sofalcone in the solvent of choice, which should be purged with an inert gas. Sofalcone is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of sofalcone in these solvents is approximately 16 and 12 mg/ml, respectively.

Sofalcone is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, sofalcone should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Sofalcone has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Sofalcone is an anti-ulcer chalcone derivative.^{1,2} It blocks inflammation by suppressing the production of nitric oxide, TNF- α , and MCP-1 while inducing the expression of heme oxygenase-1 when used at 10-20 μ M.³ Sofalcone (20-50 μ M) induces VEGF expression in gastric epithelial cells and promotes gastric ulcer healing following eradication therapy for *H. pylori*.⁴⁻⁶

References

1. Kimura, M., Saziki, R., Arai, I., et al. *Japan. J. Pharmacol.* **35**, 389-396 (1984).
2. Suwa, T., Nakazima, M., Shinozaki, A., et al. *Japan. J. Pharmacol.* **35**, 47-53 (1984).
3. Tanaka, H., Nakamura, S., Onda, K., et al. *Biochem. Biophys. Res. Commun.* **381(4)**, 566-571 (2009).
4. Isomoto, H., Furusu, H., Ohnita, K., et al. *World J. Gastroenterol.* **11(11)**, 1629-1633 (2005).
5. Higuchi, K., Watanabe, T., Tanigawa, T., et al. *J. Gastroenterol. Hepatol.* **25**, S155-S160 (2010).
6. Shibuya, A., Onda, K., Kawahara, H., et al. *Biochem. Biophys. Res. Commun.* **398(3)**, 581-584 (2010).

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, 11/07/2022

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