

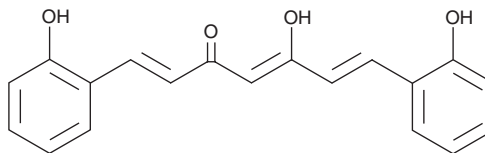
PRODUCT INFORMATION



Salicylcurcumin

Item No. 11878

CAS Registry No.: 1236545-54-2
Formal Name: (1E,4Z,6E)-5-hydroxy-1,7-bis(2-hydroxyphenyl)-1,4,6-heptatrien-3-one
Synonym: Salicylcurcuminoid
MF: C₁₉H₁₆O₄
FW: 308.3
Purity: ≥95%
UV/Vis.: λ_{max}: 288, 410 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

Salicylcurcumin is supplied as a crystalline solid. A stock solution may be made by dissolving the salicylcurcumin in the solvent of choice, which should be purged with an inert gas. Salicylcurcumin is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of salicylcurcumin in ethanol and DMSO is approximately 20 mg/ml and approximately 30 mg/ml in DMF.

Description

Curcuminoids are natural and synthetic analogs of curcumin (Item No. 81025), a natural polyphenol with modulating effects in inflammation, cancer, and immunity.¹ Salicylcurcumin is a synthetic curcuminoid which, like bisdemethoxycurcumin (Item No. 10960), lacks methoxy groups on both phenols. Furthermore, the hydroxyl groups on each phenol are positioned ortho, as in salicylates, rather than in the para position typical of curcuminoids. Salicylcurcumin has antioxidant and anticarcinogenic properties.^{2,3} Also, it increases the activity of quinone reductase in mouse hepatoma cells in a dose-dependent fashion, doubling activity at 0.3 μM salicylcurcumin.^{4,5}

References

1. Gryniewicz, G. and Slifirski, P. Curcumin and curcuminoids in quest for medicinal status. *Acta Biochim. Pol.* **59(2)**, 201-212 (2012).
2. Manju, M., Sherin, T.G., Rajasekharan, K.N., *et al.* Curcumin analogue inhibits lipid peroxidation in a freshwater teleost, *Anabas testudineus* (Bloch)--an in vitro and in vivo study. *Fish Physiol. Biochem.* **35(3)**, 413-420 (2009).
3. Anto, R.J., George, J., Babu, K.V., *et al.* Antimutagenic and anticarcinogenic activity of natural and synthetic curcuminoids. *Mutat. Res.* **370(2)**, 127-131 (1996).
4. Dinkova-Kostova, A.T. and Talalay, P. Relation of structure of curcumin analogs to their potencies as inducers of Phase 2 detoxification enzymes. *Carcinogenesis* **20(5)**, 911-914 (1999).
5. Magesh, S., Chen, Y., and Hu, L. Small molecule modulators of Keap1-Nrf2-ARE pathway as potential preventive and therapeutic agents. *Med. Res. Rev.* **32(4)**, 687-726 (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, 02/26/2024

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