PRODUCT INFORMATION



Zerumbone

Item No. 15400

CAS Registry No.:	471-05-6
Formal Name:	(2E,6E)-2,6,9,9-tetramethyl-2,6,10E-
	cycloundecatrien-1-one
MF:	$C_{15}H_{22}O$
FW:	218.3
Purity:	≥95%
Supplied as:	Solid white crystals or powder
Storage:	-20°C
Stability:	≥4 years
Item Origin:	Plant/Curcuma zerumbet
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.	



Laboratory Procedures

Zerumbone is supplied as solid white crystals or powder. A stock solution may be made by dissolving the zerumbone in the solvent of choice. Zerumbone is soluble in organic solvents such as ethanol and DMSO, which should be purged with an inert gas.

Description

Zerumbone is a natural monocyclic sesquiterpene first isolated from rhizomes of the wild ginger Z. zerumbet.¹ It potently inhibits the activation of Epstein-Barr virus by phorbol esters ($IC_{50} = 140 \text{ nM}$).^{1,2} Zerumbone inhibits Sonic hedgehog signaling and induces apoptosis in cancer cell lines.³⁻⁵ It has also been shown to have anti-oxidant and anti-inflammatory activities, contributing to immunomodulatory and hepatoprotective effects.⁶

References

- 1. Murakami, A., Takahashi, M., Jiwajinda, S., et al. Identification of zerumbone in Zingiber zerumbet smith as a potent inhibitor of 12-O-tetradecanoylphorbol-13-acetate-induced Epstein-Barr virus activation. Biosci. Biotechnol. Biochem. 63(10), 1811-1812 (1999).
- 2. Murakami, A., Takahashi, D., Kinoshita, T., et al. Zerumbone, a southeast asian ginger sesquiterpene, markedly suppresses free radical generation, proinflammatory protein production, and cancer cell proliferation accompanied by apoptosis: The α , β -unsaturated carbonyl group is a prerequisite. Carcinogenesis 23(5), 795-802 (2002).
- 3. Stanton, B.Z. and Peng, L.F. Small-molecule modulators of the Sonic Hedgehog signaling pathway. Mol. Biosyst. 6(1), 44-54 (2010).
- 4. Sun, Y., Sheng, Q., Cheng, Y., et al. Zerumbone induces apoptosis in human renal cell carcinoma via Gli-1/Bcl-2 pathway. Pharmazie 68(2), 141-145 (2013).
- 5. Sehrawat, A., Arlotti, J.A., Murakami, A., et al. Zerumbone causes Bax- and Bak-mediated apoptosis in human breast cancer cells and inhibits orthotopic xenograft growth in vivo. Breast Cancer Res. Treat. 136(2), 429-441 (2012).
- 6. Rahman, H.S., Rasedee, A., How, C.W., et al. Zerumbone-loaded nanostructured lipid carriers: Preparation, characterization, and antileukemic effect. Int. J. Nanomedicine 8, 2769-2781 (2013).

WARNING THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFFTY 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, 11/16/2022

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