

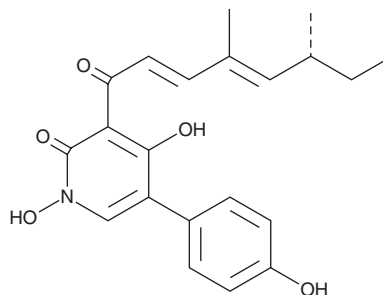
# PRODUCT INFORMATION



## Tenellin

Item No. 28090

CAS Registry No.: 53823-15-7  
Formal Name: 3-[(2E,4E,6R)-4,6-dimethyl-1-oxo-2,4-octadien-1-yl]-1,4-dihydroxy-5-(4-hydroxyphenyl)-2(1H)-pyridinone  
MF:  $C_{21}H_{23}NO_5$   
FW: 369.4  
Purity:  $\geq 95\%$   
Supplied as: A solid  
Storage:  $-20^\circ\text{C}$   
Stability:  $\geq 4$  years  
Item Origin: Fungus/*Beauveria* sp.



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

### Laboratory Procedures

Tenellin is supplied as a solid. A stock solution may be made by dissolving the tenellin in the solvent of choice. Tenellin is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide.

### Description

Tenellin is a fungal metabolite that has been found in *Beauveria*.<sup>1</sup> It inhibits  $Mg^{2+}$ -,  $Ca^{2+}$ -, and  $Na^+/K^+$ -ATPase activities in equine erythrocytes by 51, 57, and 74%, respectively, when used at a concentration of 200  $\mu\text{g}/\text{ml}$ .<sup>1</sup> Tenellin is cytotoxic to Sf9 and Sf21 insect cells with 50% cytotoxic concentration ( $CC_{50}$ ) values of 4.84 and 11.95  $\mu\text{M}$ , respectively.<sup>2</sup>

### References

1. Jeffs, L.B. and Khachatourians, G.G. Toxic properties of *Beauveria* pigments on erythrocyte membranes. *Toxicon*. **35(8)**, 1351-1356 (1997).
2. Arboleda Valencia, J.W., Gaitán Bustamante, A.L., Jiménez, A.V., *et al.* Cytotoxic activity of fungal metabolites from the pathogenic fungus *Beauveria bassiana*: An intraspecific evaluation of beauvericin production. *Curr. Microbiol.* **63(3)**, 306-312 (2011).

#### 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, 08/11/2023

#### 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