

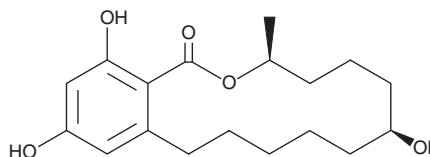
PRODUCT INFORMATION



Taleranol

Item No. 28299

CAS Registry No.: 42422-68-4
Formal Name: (3S,7S)-3,4,5,6,7,8,9,10,11,12-decahydro-7,14,16-trihydroxy-3-methyl-1H-2-benzoxacyclotetradecin-1-one
Synonyms: P 1560, β -Zearalanol, β -Zearanol
MF: C₁₈H₂₆O₅
FW: 322.4
Purity: \geq 98%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

Taleranol is supplied as a solid. A stock solution may be made by dissolving the taleranol in the solvent of choice, which should be purged with an inert gas. Taleranol is soluble in the organic solvent methanol at a concentration of approximately 10 mg/ml.

Description

Taleranol is a β -resorcylic acid lactone and an isomer of α -zearalanol (Item No. 23424).¹ *In vivo*, taleranol (1-500 μ g/animal) induces synthesis of the uterine protein IP and increases uterine weight in rats.

Reference

1. Katzenellenbogen, B.S., Katzenellenbogen, J.A., and Mordecai, D. Zearalenones: Characterization of the estrogenic potencies and receptor interactions of a series of fungal β -resorcylic acid lactones. *Endocrinology* **105**(1), 33-40 (1979).

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, 10/04/2022

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