

# PRODUCT INFORMATION

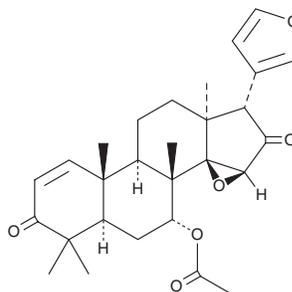


## Epoxyzadiradione

Item No. 44540

**CAS Registry No.:** 18385-59-6  
**Formal Name:** (5 $\alpha$ ,7 $\alpha$ ,13 $\alpha$ ,14 $\beta$ ,15 $\beta$ ,17 $\alpha$ )-7-(acetyloxy)-14,15:21,23-diepoxy-4,4,8-trimethyl-24-norchola-1,20,22-triene-3,16-dione

**MF:** C<sub>28</sub>H<sub>34</sub>O<sub>6</sub>  
**FW:** 466.6  
**Purity:**  $\geq$ 98%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:**  $\geq$ 4 years  
**Item Origin:** Synthetic



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

### Laboratory Procedures

Epoxyzadiradione is supplied as a solid. A stock solution may be made by dissolving the epoxyzadiradione in the solvent of choice, which should be purged with an inert gas. Epoxyzadiradione is slightly soluble (0.1-1 mg/ml) in acetonitrile and chloroform.

### Description

Epoxyzadiradione is a limonoid that has been found in *A. indica* and has diverse biological activities.<sup>1,2</sup> It inhibits macrophage migration inhibitor factor (MIF) tautomerase activity (IC<sub>50</sub>s = 6.4, 9.5, and 9.8  $\mu$ M for the human, *P. falciparum*, and *P. yoelii* enzymes, respectively).<sup>1</sup> Epoxyzadiradione (400  $\mu$ M) inhibits tau aggregation and induces disaggregation of preformed tau aggregates in cell-free assays.<sup>1</sup> *In vivo*, epoxyzadiradione (25, 50, and 100 mg/kg) prevents LPS- and *P. yoelii* MIF-induced IL-6, IL-1 $\alpha$ , IL-1 $\beta$ , and TNF- $\alpha$  production in mice.<sup>1</sup>

### References

1. Alam, A., Haldar, S., Thulasiram, H.V., *et al.* Novel anti-inflammatory activity of epoxyzadiradione against macrophage migration inhibitory factor: Inhibition of tautomerase and proinflammatory activities of macrophage migration inhibitory factor. *J. Biol. Chem.* **287**(29), 24844-24861 (2012).
2. Gorantla, N.V., Das, R., Chidambaram, H., *et al.* Basic limonoid modulates chaperone-mediated proteostasis and dissolve Tau fibrils. *Sci. Rep.* **10**(1), 4023 (2020).

#### 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, 12/17/2025

#### 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