

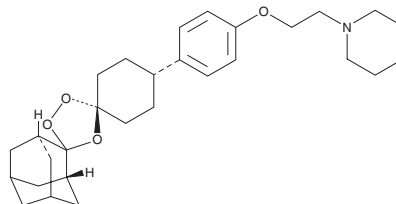
# PRODUCT INFORMATION



## Artefenomel

Item No. 30677

**CAS Registry No.:** 1029939-86-3  
**Formal Name:** 4-[2-(4-*cis*-dispiro[cyclohexane-1,3'-[1,2,4]trioxolane-5',2''-tricyclo[3.3.1.1<sup>3,7</sup>]decan]-4-ylphenoxy)ethyl]-morpholine  
**Synonym:** OZ439  
**MF:** C<sub>28</sub>H<sub>39</sub>NO<sub>5</sub>  
**FW:** 469.6  
**Purity:** ≥98%  
**Supplied as:** A 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

Artefenomel is supplied as a solid. A stock solution may be made by dissolving the artefenomel in the solvent of choice, which should be purged with an inert gas. Artefenomel is soluble in the organic solvent DMSO at a concentration of approximately 10 mM.

### Description

Artefenomel is an ozonide antimalarial agent.<sup>1</sup> It is active against 3D7, Cam3.II, and Cam3.II\_rev *P. falciparum* strains (LC<sub>50</sub>s = 8.7, 4.4, and 6.1 nM, respectively). Artefenomel reduces the number of trophozoites in *P. falciparum*-infected isolated human red blood cells (IC<sub>50</sub> = 31 nM).<sup>2</sup> *In vivo*, artefenomel (30 mg/kg) is 100% curative in a mouse model of *P. berghei* infection.<sup>3</sup>

### References

1. Yang, T., Xie, S.C., Cao, P., *et al.* Comparison of the exposure time dependence of the activities of synthetic ozonide antimalarials and dihydroartemisinin against K13 wild-type and mutant *Plasmodium falciparum* strains. *Antimicrob. Agents Chemother.* **60(8)**, 4501-4510 (2016).
2. Giannangelo, C., Stingelin, L., Yang, T., *et al.* Parasite-mediated degradation of synthetic ozonide antimalarials impacts *in vitro* antimalarial activity. *Antimicrob. Agents Chemother.* **62(3)**, e01566-e01617 (2018).
3. Kim, H.S., Hammill, J.T., and Guy, R.K. Seeking the elusive long-acting ozonide: Discovery of artefenomel (OZ439). *J. Med. Chem.* **60(7)**, 2651-2653 (2017).

#### 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, 10/14/2022

#### 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