

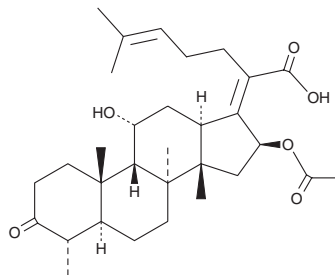
PRODUCT INFORMATION



3-keto Fusidic Acid

Item No. 33855

CAS Registry No.: 4680-37-9
Formal Name: (4 α ,8 α ,9 β ,11 α ,13 α ,14 β ,16 β ,17Z)-16-(acetyloxy)-11-hydroxy-3-oxo-29-nordammara-17(20),24-dien-21-oic acid
Synonyms: 3-keto FA, 3-Oxofusidic Acid
MF: C₃₁H₄₆O₆
FW: 514.7
Purity: \geq 95%
Supplied as: A solid
Storage: -20°C
Stability: \geq 2 years
Item Origin: Synthetic



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

Laboratory Procedures

3-keto Fusidic acid (3-keto FA) is supplied as a solid. A stock solution may be made by dissolving the 3-keto FA in the solvent of choice, which should be purged with an inert gas. 3-keto FA is slightly soluble in chloroform and methanol.

Description

3-keto FA is an active metabolite of the antibiotic fusidic acid (Item No. 14825).¹ 3-keto FA is active against *M. tuberculosis* (MIC₉₉ = 1.25 μ M).

Reference

1. Strydom, N., Kaur, G., Dziwornu, G.A., *et al.* Pharmacokinetics and organ distribution of C 3 alkyl esters as potential antimycobacterial prodrugs of fusidic acid. *ACS Infect. Dis.* **6(3)**, 459-466 (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, 04/07/2021

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