

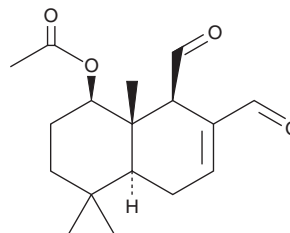
PRODUCT INFORMATION



1 β -Acetoxypolygodial

Item No. 29901

CAS Registry No.: 1818425-44-3
Formal Name: (1R,4aS,8R,8aS)-8-(acetyloxy)-
1,4,4a,5,6,7,8,8a-octahydro-5,5,8a-trimethyl-
1,2-naphthalenedicarboxaldehyde
MF: C₁₇H₂₄O₄
FW: 292.4
Purity: $\geq 95\%$
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥ 4 years
Item Origin: Plant/*Tasmannia lanceolata*



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

Laboratory Procedures

1 β -Acetoxypolygodial is supplied as a crystalline solid. A stock solution may be made by dissolving the 1 β -acetoxypolygodial in the solvent of choice, which should be purged with an inert gas. 1 β -Acetoxypolygodial is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide.

Description

1 β -Acetoxypolygodial is a sesquiterpenoid that has been found in *T. lanceolata* and has antiprotozoal activity.^{1,2} It is active against *T. cruzi* amastigotes and trypomastigotes (GI₅₀s = 57.8 and 72 μ M, respectively) but not *T. cruzi* epimastigotes (GI₅₀ = >100 μ M).²

References

1. Deans, R.M., Gardiner, M.G., Horne, J., *et al.* Isolation and characterization of 1 β -acetoxypolygodial from *Tasmannia lanceolata*. *Asian J. Chem.* **3**(11), 1193-1196 (2014).
2. Turner, D.N., Just, J., Dasari, R., *et al.* Activity of natural and synthetic polygodial derivatives against *Trypanosoma cruzi* amastigotes, trypomastigotes and epimastigotes. *Nat. Prod. Res.* (2019).

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, 10/21/2022

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