

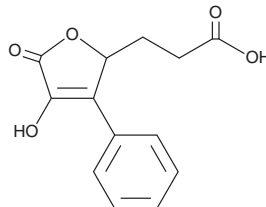
PRODUCT INFORMATION



WF-3681

Item No. 33523

CAS Registry No.: 105364-56-5
Formal Name: 2,5-dihydro-4-hydroxy-5-oxo-3-phenyl-furanpropanoic acid
Synonym: (±)-WF-3681
MF: C₁₃H₁₂O₅
FW: 248.2
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Fungus/*Chaetomella* sp.



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

Laboratory Procedures

WF-3681 is supplied as a solid. A stock solution may be made by dissolving the WF-3681 in the solvent of choice, which should be purged with an inert gas. WF-3681 is soluble in methanol and DMSO.

Description

WF-3681 is a butenolide fungal metabolite that has been found in *Acremonium*.¹ It inhibits aldose reductase (IC₅₀ = 0.25 μM for the rabbit lens enzyme).²

References

1. Ghisalberti, E.L., Hargreaves, J.R., and McConville, E. Butenolides from a cultured microfungus, *Acremonium* sp. *Nat. Prod. Res.* **18(2)**, 105-109 (2004).
2. Nishikawa, M., Yoshida, K., Okamoto, M., et al. Studies on WE-3681, a novel aldose reductase inhibitor. III. Effects of WF-3681 and its derivatives on sorbitol accumulation in diabetic rats. *J. Antibiot. (Tokyo)* **43(9)**, 1186-1188 (1990).

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, 08/25/2023

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