

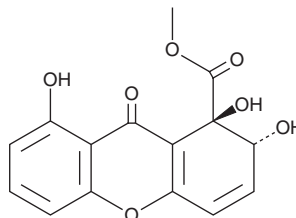
PRODUCT INFORMATION



Globosuxanthone A

Item No. 27486

CAS Registry No.: 917091-74-8
Formal Name: (1R,2R)-2,9-dihydro-1,2,8-trihydroxy-9-oxo-1H-xanthene-1-carboxylic acid, methyl ester
MF: $C_{15}H_{12}O_7$
FW: 304.3
Purity: $\geq 95\%$
Supplied as: A solid
Storage: -20°C
Stability: ≥ 4 years
Item Origin: Fungus/Unidentified sp.



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

Laboratory Procedures

Globosuxanthone A is supplied as a solid. A stock solution may be made by dissolving the globosuxanthone A in the solvent of choice. Globosuxanthone A is soluble in organic solvents such as ethanol, methanol, and DMSO, which should be purged with an inert gas. The solubility of globosuxanthone A in ethanol is approximately 1 mg/ml.

Description

Globosuxanthone A is a fungal metabolite originally isolated from *C. globosum*.¹ It is cytotoxic to various human solid tumor cell lines, including MCF-7, PC3, LNCap, and DU145 cells (IC_{50} s = 1.3, 0.65, 1.5, and 1.2 μM , respectively). It induces accumulation of cells in the G_2/M and S phases of the cell cycle in NCI-H460 and PC3M cells. Globosuxanthone A is also active against *C. albicans* but not *S. aureus*, *E. coli*, or *M. hiemalis*.²

References

1. Wijeratne, E.M.K., Turbyville, T.J., Fritz, A., *et al.* A new dihydroxanthone from a plant-associated strain of the fungus *Chaetomium globosum* demonstrates anticancer activity. *Bioorg. Med. Chem.* **14**(23), 7917-7923 (2006).
2. Yamazaki, H., Rotinsulu, H., Kaneko, T., *et al.* A new dibenz[b,e]oxepine derivative, 1-hydroxy-10-methoxy-dibenz[b,e]oxepin-6,11-dione, from a marine-derived fungus, *Beauveria bassiana* TPU942. *Mar. Drugs* **10**(12), 2691-2697 (2012).

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, 08/09/2023

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