

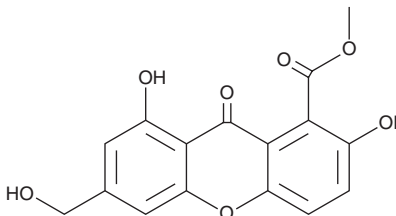
# PRODUCT INFORMATION



## Sydowinin B

Item No. 31670

**CAS Registry No.:** 58450-00-3  
**Formal Name:** 2,8-dihydroxy-6-(hydroxymethyl)-9-oxo-9H-xanthene-1-carboxylic acid, methyl ester  
**Synonym:** MS-347b  
**MF:** C<sub>16</sub>H<sub>12</sub>O<sub>7</sub>  
**FW:** 316.3  
**Purity:** ≥95%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years  
**Item Origin:** Fungus/*Aspergillus sydowii*



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

### Laboratory Procedures

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

### Description

Sydowinin B is a xanthone polyketide that has been found in *A. sydowii* and has immunosuppressant activity.<sup>1</sup> It inhibits LPS- or concanavalin A-induced proliferation of isolated mouse splenic lymphocytes (IC<sub>50</sub>s = 19.2 and 20.8 µg/ml, respectively).<sup>2</sup> Sydowinin B inhibits superoxide generation in, and elastase release from, isolated human neutrophils induced by cytochalasin B (Item No. 11328) and N-Formyl-Met-Leu-Phe (fMLF; Item No. 21495) with IC<sub>50</sub> values of 21.2 and 12.62 µM, respectively.<sup>3</sup>

### References

1. Hamasaki, T., Sato, Y., and Hatsuda, Y. Structure of sydowinin A, sydowinin B, and sydowinol, metabolites from *Aspergillus sydowii*. *Agr. Biol. Chem.* **39**(12), 2341-2345 (1975).
2. Liu, H., Chen, S., Liu, W., *et al.* Polyketides with immunosuppressive activities from mangrove endophytic fungus *Penicillium* sp. ZJ-SY<sub>2</sub>. *Mar. Drugs* **14**(12), 217 (2016).
3. Chung, Y.-M., Wei, C.-K., Chuang, D.-W., *et al.* An epigenetic modifier enhances the production of anti-diabetic and anti-inflammatory sesquiterpenoids from *Aspergillus sydowii*. *Bioorg. Med. Chem.* **21**(13), 3866-3872 (2013).

#### 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, 08/11/2023

#### 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](http://WWW.CAYMANCHEM.COM)