

# PRODUCT INFORMATION



## Stachybotrysin B

Item No. 29583

**CAS Registry No.:** 2098376-42-0  
**Formal Name:** (1'R,2'R,4'aS,6'R,8'aS)-6'-(acetyloxy)-3',4',4'a,5',6',7',8',8'a-octahydro-4-hydroxy-6-(hydroxymethyl)-2',5',5',8'a-tetramethyl-spiro[benzofuran-2(3H),1'(2'H)-naphthalene]-7-carboxaldehyde

**MF:** C<sub>25</sub>H<sub>34</sub>O<sub>6</sub>

**FW:** 430.5

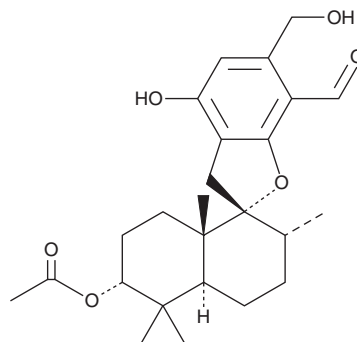
**Purity:** ≥95%

**Supplied as:** A solid

**Storage:** -20°C

**Stability:** ≥4 years

**Item Origin:** Fungus/*Stachybotrys* sp.



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

### Laboratory Procedures

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

### Description

Stachybotrysin B is a fungal metabolite originally isolated from *S. chartarum* and has antiviral and anticancer activities.<sup>1,2</sup> It has antiviral activity against HIV in SupT1 cells (IC<sub>50</sub> = 19.2 μM).<sup>1</sup> Stachybotrysin B is cytotoxic to K562, HeLa, and HL-60 cells (IC<sub>50</sub>s = 21.72, 39.63, and 18.5 μM, respectively).<sup>2</sup>

### References

1. Zhao, J., Feng, J., Tan, Z., *et al.* Stachybotrysin A-G, phenylspirodrimane derivatives from the fungus *Stachybotrys chartarum*. *J. Nat. Prod.* **80**(6), 1819-1826 (2017).
2. Ma, X.-h., Zheng, W.-m., Sun, K.-h., *et al.* Two new phenylspirodrimanes from the deep-sea derived fungus *Stachybotrys* sp. MCCC 3A00409. *Nat. Prod. Res.* **33**(3), 386-392 (2018).

#### 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, 08/15/2023

#### 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)