

# PRODUCT INFORMATION



## Stachybotrylactam

Item No. 22584

CAS Registry No.: 163391-76-2

Formal Name: (1'R,2'R,4'aS,6'R,8'aS)-3',4',4'a,5',6',7,7',8,8'a-decahydro-4,6'-dihydroxy-2',5',5',8'a-tetramethyl-spiro[2H-furo[2,3-e]isoindole-2,1'(2'H)-naphthalen]-6(3H)-one

MF: C<sub>23</sub>H<sub>31</sub>NO<sub>4</sub>

FW: 385.5

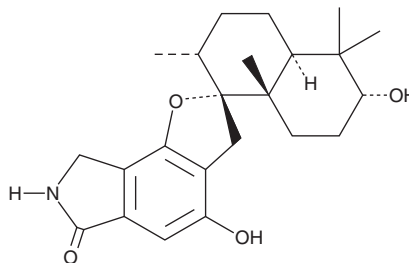
Purity: ≥95%

UV/Vis.: λ<sub>max</sub>: 219 nm

Supplied as: A lyophilisate

Storage: -20°C

Stability: ≥4 years



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

### Laboratory Procedures

Stachybotrylactam is supplied as a lyophilisate. A stock solution may be made by dissolving the stachybotrylactam in the solvent of choice, which should be purged with an inert gas. Stachybotrylactam is soluble in ethanol, methanol, DMSO, and dimethyl formamide.

### Description

Stachybotrylactam is a spirodihydrobenzofuran lactam mycotoxin originally isolated from *Stachybotrys* with antiviral activity.<sup>1,2</sup> It inhibits HIV-1 protease *in vitro* with an IC<sub>50</sub> value of 161 μM.<sup>2</sup> Stachybotrylactam has no effect on oleic acid-elicited intracellular lipid accumulation and has no observable cytotoxicity at concentrations up to 100 μM in HepG2 cells.<sup>3</sup>

### References

1. Jarvis, B.B., Salemme, J., Morais, A., *et al.* *Stachybotrys* toxins. 1. *Nat. Toxins.* **3**(1), 10-16 (1995).
2. Roggo, B.E., Petersen, F., Sills, M., *et al.* Novel spirodihydrobenzofuran lactams as antagonists of endothelin and as inhibitors of HIV-1 protease produced by *Stachybotrys* Sp. I. Fermentation, isolation and biological activity. *J. Antibiot. (Tokyo)* **49**(1), 13-19 (1996).
3. Li, Y., Wu, C., Liu, D., *et al.* Chartaractams A-P, phenylspirodrimanones from the sponge-associated fungus *Stachybotrys chartarum* with antihyperlipidemic activities. *J. Nat. Prod.* **77**(1), 138-147 (2014).

#### 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, 11/11/2022

#### 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