

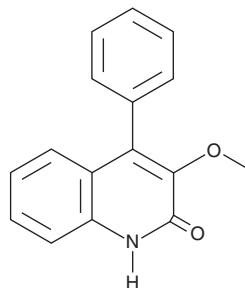
# PRODUCT INFORMATION



## O-Methylviridicatin

Item No. 19446

**CAS Registry No.:** 6152-57-4  
**Formal Name:** 3-methoxy-4-phenyl-2(1H)-quinolinone  
**Synonyms:** AIDS-089094, 3-methoxy-4-phenyl-Carbostyryl, Viridicatin methyl ester  
**MF:** C<sub>16</sub>H<sub>13</sub>NO<sub>2</sub>  
**FW:** 251.3  
**Purity:** ≥99%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years  
**Special Conditions:** XXXX



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

### Laboratory Procedures

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

O-Methylviridicatin is sparingly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

### Description

O-Methylviridicatin is a natural derivative of the alkaloid mycotoxin viridicatin, which is produced by several species of *Penicillium*.<sup>1</sup> O-Methylviridicatin blocks activation of the HIV long terminal repeat by TNF- $\alpha$  in HeLa cells (IC<sub>50</sub> = 5  $\mu$ M).<sup>2</sup> It also inhibits virus production in OM-10.1 cells (HL-60 promyelocytes infected with HIV-1) treated with TNF- $\alpha$  (IC<sub>50</sub> = 2.5  $\mu$ M).<sup>2</sup>

### References

1. Santini, A., Mikusovj, P., Sulyok, M., *et al.* *Penicillium* strains isolated from Slovak grape berries taxonomy assessment by secondary metabolite profile. *Mycotoxin Res.* **30**(4), 213-220 (2014).
2. Heguy, A., Cai, P., Meyn, P., *et al.* Isolation and characterization of the fungal metabolite 3-O-methylviridicatin as an inhibitor of tumour necrosis factor  $\alpha$ -induced human immunodeficiency virus replication. *Antivir. Chem. Chemother.* **9**(2), 149-155 (1998).

#### 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, 12/15/2022

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