

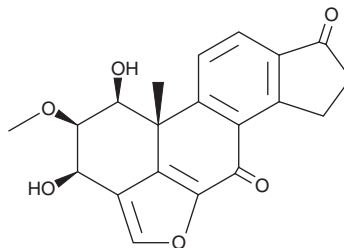
# PRODUCT INFORMATION



## Viridiol

Item No. 19607

CAS Registry No.: 23820-80-6  
Formal Name: 1 $\beta$ ,3 $\beta$ -dihydroxy-2 $\beta$ -methoxy-  
18-norandrosta-5,8,11,13-  
tetraeno[6,5,4-bc]furan-7,17-dione  
Synonym: NSC 658706  
MF: C<sub>20</sub>H<sub>18</sub>O<sub>6</sub>  
FW: 354.4  
Purity:  $\geq 95\%$   
Supplied as: A solid  
Storage: -20°C  
Stability:  $\geq 4$  years  
Special Conditions: Keep cool and dry.  
Item Origin: Fungus/*Trichoderma* sp.



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

### Laboratory Procedures

Viridiol is supplied as a solid. A stock solution may be made by dissolving the viridiol in the solvent of choice. Viridiol is soluble in the organic solvent DMSO.

### Description

Viridiol is a fungal steroid that demonstrates phytotoxic activity.<sup>1</sup> It is a metabolite of the fungistatic compound, viridin. Both viridin and viridiol are reported to inhibit PI3K.<sup>2</sup>

### References

1. Andersson, P.F., Johansson, S.B.K., Stenlid, J., *et al.* Isolation, identification and necrotic activity of viridiol from *Chalara fraxinea*, the fungus responsible for dieback of ash. *For. Path.* **40**(1), 43-46 (2009).
2. Dodge, J.A., Sato, M., and Vlahos, C.J. Inhibition of phosphatidylinositol 3-kinase with viridin, demethoxyviridin, viridiol, demethoxyviridiol, virone, wortmannolone, and analogs thereof. 5,726,167 (1998), 805,704.

#### 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