

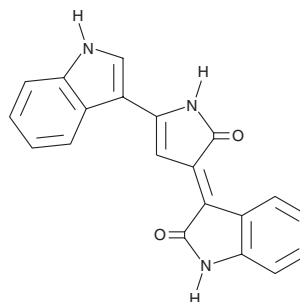
PRODUCT INFORMATION



Deoxyviolacein

Item No. 23477

CAS Registry No.: 5839-61-2
Formal Name: (3E)-3-[1,2-dihydro-5-(1H-indol-3-yl)-2-oxo-3H-pyrrol-3-ylidene]-1,3-dihydro-2H-indol-2-one
MF: C₂₀H₁₃N₃O₂
FW: 327.3
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

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

Deoxyviolacein 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

Deoxyviolacein is a bacterial metabolite and byproduct in the biosynthesis of the bisindole alkaloid violacein (Item No. 27959) that has anticancer, antibacterial, and antifungal properties.^{1,2} It inhibits proliferation of hepatocellular carcinoma cells when used at concentrations ranging from 0.1 to 1 μM.² Deoxyviolacein (125 μg/ml) has antibacterial activity against Gram-positive bacteria, including *S. aureus*, *B. subtilis*, and *B. megaterium*.¹ It also has antifungal activity against *R. solani* when used at a concentration of 2 mg/ml.

References

1. Wang, H., Wang, F., Zhu, X., *et al.* Biosynthesis and characterization of violacein, deoxyviolacein and oxyviolacein in heterologous host, and their antimicrobial activities. *Biochem. Eng. J.* **67**, 148-155 (2012).
2. Jiang, P.-X., Wang, H.-S., Xiao, S., *et al.* Pathway redesign for deoxyviolacein biosynthesis in *Citrobacter freundii* and characterization of this pigment. *Appl. Microbiol. Biotechnol.* **94**(6), 1521-1532 (2012).

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, 02/19/2025

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