

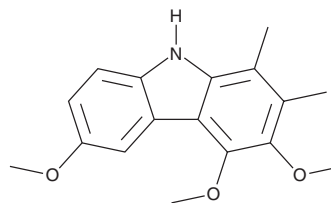
PRODUCT INFORMATION



Carbazomycin D

Item No. 33091

CAS Registry No.: 108073-63-8
Formal Name: 3,4,6-trimethoxy-1,2-dimethyl-9H-carbazole
MF: C₁₇H₁₉NO₃
FW: 285.3
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Bacterium/*Streptomyces* sp.



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

Laboratory Procedures

Carbazomycin D is supplied as a solid. A stock solution may be made by dissolving the carbazomycin D in the solvent of choice, which should be purged with an inert gas. Carbazomycin D is soluble in methanol and DMSO.

Description

Carbazomycin D is a bacterial metabolite that has been found in *Streptomyces* and has diverse biological activities.^{1,2} It is active against the fungi *T. asteroides* and *T. mentagrophytes* (MIC = 100 µg/ml for both) and the bacterium *M. tuberculosis* (IC₅₀ = 25 µg/ml). Carbazomycin D is cytotoxic to MCF-7, KB, NCI H187, and Vero cells (IC₅₀s = 21.3, 33.2, 12.9, and 34.3 µg/ml, respectively).²

References

1. Naid, T., Kitahara, T., Kaneda, M., *et al.* Carbazomycins C, D, E and F, minor components of the carbazomycin complex. *J. Antibiot. (Tokyo)* **40(2)**, 157-164 (1987).
2. Intaraudom, C., Rachtawee, P., Suvannakad, R., *et al.* Antimalarial and antituberculosis substances from *Streptomyces* sp. BCC26924. *Tetrahedron* **67(39)**, 7593-7597 (2011).

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