

PRODUCT INFORMATION



SCH 38519

Item No. 28092

CAS Registry No.: 114970-20-6

Formal Name: [3aR-(3a α ,5 α ,9 β ,10 β ,11 α ,13 β ,14b α)]-11-(dimethylamino)-3,3a,5,10,11,12,13,14b-octahydro-7,10-dihydroxy-5,9-dimethyl-9,13-epoxycyclohepta[7,8]naphtho[2,3-d]furo[3,2-b]pyran-2,6,14(9H)-trione

MF: C₂₄H₂₅NO₈

FW: 455.5

Purity: $\geq 85\%$

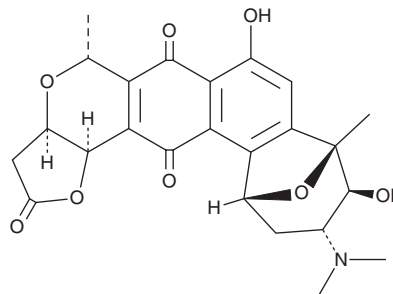
UV/Vis.: $\lambda_{\max}$: 218 nm

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

SCH 38519 is supplied as a solid. A stock solution may be made by dissolving the SCH 38519 in the solvent of choice, which should be purged with an inert gas. SCH 38519 is soluble in organic solvents such as ethanol, methanol, DMSO, and dimethyl formamide.

Description

SCH 38519 is an isochromanequinone fungal metabolite originally isolated from *Thermomonospora*.¹ It inhibits aggregation induced by thrombin (Item No. 13188), as well as inhibits serotonin secretion in human platelets (IC₅₀s = 68 and 61 $\mu\text{g/ml}$, respectively).² It inhibits the growth of various Gram-negative and Gram-positive bacteria (mean MICs = 122.1 and 0.92 $\mu\text{g/ml}$, respectively).

References

1. Hegde, V.R., King, A.H., Patel, M.G., *et al.* A new isochromanequinone antibiotic, SCH 38519, with a novel sugar linkage. *Tetrahedron Lett.* **28**(39), 4485-4488 (1987).
2. Patel, M., Hegde, V., Horan, A., *et al.* Sch 38519, a novel platelet aggregation inhibitor produced by a *Thermomonospora* sp. Taxonomy, fermentation, isolation, physico-chemical properties, structure and biological properties. *J. Antibiot. (Tokyo)* **42**(7), 1063-1069 (1989).

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, 07/22/2025

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