

PRODUCT INFORMATION

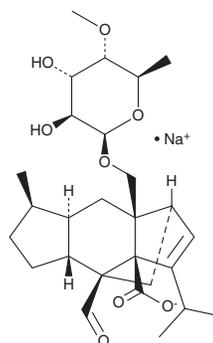


Sordarin (sodium salt)

Item No. 26255

CAS Registry No.: 463356-00-5
Formal Name: (1R,3aR,4S,4aR,7R,7aR,8aS)-8a-[[[(6-deoxy-4-O-methyl-β-D-altropyranosyl)oxy]methyl]-4-formyl-4,4a,5,6,7,7a,8,8a-octahydro-7-methyl-3-(1-methylethyl)-1,4-methano-s-indacene-3a(1H)-carboxylic acid, monosodium salt

MF: C₂₇H₃₉O₈ • Na
FW: 514.6
Purity: ≥98%
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

Sordarin (sodium salt) is supplied as a solid. A stock solution may be made by dissolving the sordarin (sodium salt) in water. We do not recommend storing the aqueous solution for more than one day.

Description

Sordarin is an inhibitor of fungal protein synthesis originally isolated from *S. araneosa*.¹ It binds to elongation factor 2 (EF-2) in the presence of ribosomes and inhibits the uncoupled GTPase activity of equimolar mixtures of EF-2 and ribosomes from *C. albicans* (IC₅₀ = 0.1 μM). Sordarin inhibits protein synthesis in cell-free lysates of *C. albicans*, *C. glabrata*, and *C. neoformans* (IC₅₀s = 0.01, 0.2, and 0.06 μg/ml, respectively) but not in rabbit reticulocytes (IC₅₀ = >100 μg/ml).² It inhibits the growth of *C. albicans* (MIC = 8 μg/ml) but not *C. glabrata* or *C. neoformans* (MICs = >125 μg/ml).

References

1. Domínguez, J.M., Gómez-Lorenzo, M.G., and Martín, J.J. Sordarin inhibits fungal protein synthesis by blocking translocation differently to fusidic acid. *J. Biol. Chem.* **274**(32), 22423-22427 (1999).
2. Domínguez, J.M., Kelly, V.A., Kinsman, O.S., et al. Sordarins: A new class of antifungals with selective inhibition of the protein synthesis elongation cycle in yeasts. *Antimicrob. Agents Chemother.* **42**(9), 2274-2278 (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, 12/19/2022

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