

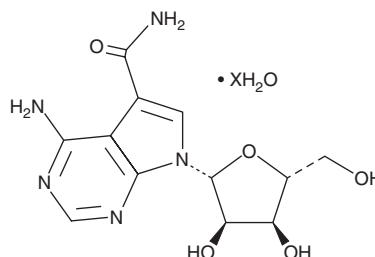
PRODUCT INFORMATION



Sangivamycin (hydrate)

Item No. 21343

Formal Name: 4-amino-7-β-D-ribofuranosyl-7H-pyrrolo[2,3-d]pyrimidine-5-carboxamide, hydrate
Synonym: NSC 65346
MF: C₁₂H₁₅N₅O₅ • XH₂O
FW: 309.3
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

Sangivamycin (hydrate) is supplied as a solid. A stock solution may be made by dissolving the sangivamycin (hydrate) in the solvent of choice, which should be purged with an inert gas. Sangivamycin (hydrate) is slightly soluble (0.1-1 mg/ml) in acetonitrile and DMSO.

Sangivamycin (hydrate) is slightly 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

Sangivamycin is a nucleoside analog.¹ It incorporates into severe acute respiratory syndrome coronavirus 2 (SARS-CoV-2) RNA and reduces the number of plaque-forming units (PFUs) in SARS-CoV-2-infected Vero E6 cells when used at a concentration of 72 nM. Sangivamycin (0.3 μM) induces cell cycle arrest at the G₂/M phase and apoptosis in MCF-7/adr cells, effects that can be prevented by the JNK inhibitor SP600125 (Item No. 10010466) or protein kinase Cδ (PKCδ) inhibitor rottlerin (Item No. 12006).²

References

1. Bennett, R.P., Yoluç, Y., Salter, J.D., *et al.* Sangivamycin is preferentially incorporated into viral RNA by the SARS-CoV-2 polymerase. *Antiviral Res.* **218**, 105716 (2023).
2. Lee, S.A. and Jung, M. The nucleoside analog sangivamycin induces apoptotic cell death in breast carcinoma MCF7/adriamycin-resistant cells via protein kinase Cdelta and JNK activation. *J. Biol. Chem.* **282**(20), 15271-15283 (2007).

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