

# PRODUCT INFORMATION



## Kirromycin

Item No. 25189

CAS Registry No.: 50935-71-2

Formal Name: (αS,2R,3R,4R,6S)-N-[(2E,4E,6S,7R)-7-[(2S,3S,4R,5R)-5-[(1E,3E,5E)-7-(1,2-dihydro-4-hydroxy-2-oxo-3-pyridinyl)-6-methyl-7-oxo-1,3,5-heptatrien-1-yl]tetrahydro-3,4-dihydroxy-2-furanyl]-6-methoxy-5-methyl-2,4-octadien-1-yl]-α-ethyltetrahydro-2,3,4-trihydroxy-5,5-dimethyl-6-(1E,3Z)-1,3-pentadien-1-yl-2H-pyran-2-acetamide

Synonym: NSC 316094

MF: C<sub>43</sub>H<sub>60</sub>N<sub>2</sub>O<sub>12</sub>

FW: 796.9

Purity: ≥95%

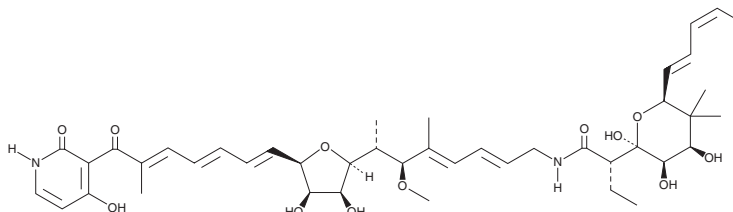
Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Item Origin: Bacterium/*Streptomyces collinus*

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



### Laboratory Procedures

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

### Description

Kirromycin is an antibiotic originally isolated from *Streptomyces* and an inhibitor of protein biosynthesis.<sup>1</sup> It inhibits isoleucine incorporation, polyphenylalanine synthesis, and growth of *B. brevis*. Kirromycin inhibits elongation factor Tu-dependent peptidyl transfer activity in *E. coli*.<sup>2</sup>

### References

1. Wolf, H., Zähler, H., and Nierhaus, K. Kirromycin, an inhibitor of the 30 S ribosomal subunits function. *FEBS Lett.* **21(3)**, 347-350 (1972).
2. Wolf, H., Chinali, G., and Parmeggiani, A. Kirromycin, an inhibitor of protein biosynthesis that acts on elongation factor Tu. *Proc. Natl. Acad. Sci. USA* **71(12)**, 4910-4914 (1974).

#### 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, 12/20/2022

#### 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