

# PRODUCT INFORMATION



## Moenomycin Complex

Item No. 15506

CAS Registry No.: 11015-37-5

Synonyms: Bambermycin, Flavomyacin, Flavophospholipol, Gainpro

MF:  $C_{69}H_{107}N_4O_{35}P$  (for A)

FW: 1,583.6 (for A)

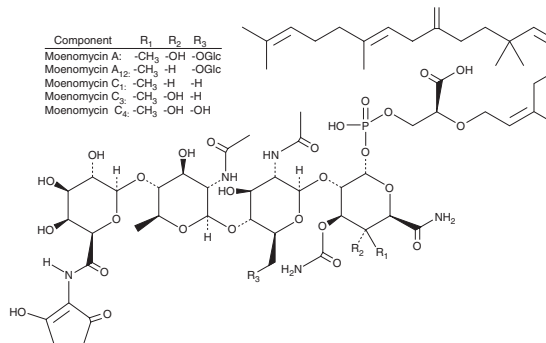
Purity:  $\geq 95\%$

Supplied as: A solid

Storage:  $-20^{\circ}C$

Stability:  $\geq 2$  years

Item Origin: Bacterium/*Streptomyces* sp.



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

### Laboratory Procedures

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

### Description

Moenomycin complex is a mixture of moenomycins A, A<sub>12</sub>, C<sub>1</sub>, C<sub>3</sub> and C<sub>4</sub>, which are antibiotics isolated from several strains of *Streptomyces* that directly inhibit bacterial peptidoglycan glycosyltransferases.<sup>1</sup> The minimal inhibitory concentration of moenomycin A against various Gram-positive bacteria ranges from 1-100 nM.<sup>1</sup>

### Reference

- Ostash, B. and Walker, S. Moenomycin family antibiotics: Chemical synthesis, biosynthesis, biological activity. *Nat. Prod. Rep.* **27(11)**, 1594-1617 (2010).

#### 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, 02/17/2021

#### 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