

# PRODUCT INFORMATION

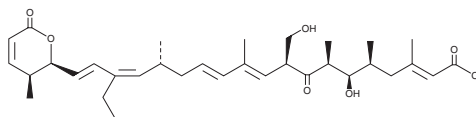


## Kazusamycin A

Item No. 20137

CAS Registry No.: 92090-94-3

Formal Name: (2E,5S,6R,7S,9S,10E,12E,15R,16Z,18E)-19-[(2S,3S)-3,6-dihydro-3-methyl-6-oxo-2H-pyran-2-yl]-17-ethyl-6-hydroxy-9-(hydroxymethyl)-3,5,7,11,15-pentamethyl-8-oxo-2,10,12,16,18-nonadecapentaenoic acid



Synonyms: CL 1,957B, Hydroxylactocin, Hydroxyleptomycin B, PD 114721

MF:  $C_{33}H_{48}O_7$

FW: 556.7

Purity:  $\geq 95\%$

Supplied as: A film

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

Stability:  $\geq 4$  years

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

### Laboratory Procedures

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

Kazusamycin A is sparingly 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

Kazusamycin A is an antibiotic from *Streptomyces* and a hydroxy analog of leptomycin B (Item No. 10004976) that demonstrates cytotoxic activity against various human and mouse tumor lines.<sup>1</sup> Additionally, at nanomolar concentrations, it inhibits nuclear export and translocation of REV, a regulatory gene product in the HIV genome.<sup>2,3</sup>

### References

1. Tunac, J. B., Graham, B. D., Dobson, W. E., *et al.* Novel antitumor antibiotics, CI-940 (PD 114,720) and PD 114,721. Taxonomy, fermentation and biological activity. *J. Antibiot. (Tokyo)* **38**(4), 460-465 (1985).
2. Wang, Y., Ponelle, M., Sanglier, J.-J., *et al.* Novel leptomycins from a *Streptomyces* strain A92-308902: Inhibitors of the nucleo-cytoplasmic translocation of the HIV-1 regulatory protein Rev. *Helv. Chim. Acta.* **80**(7), 2157-2167 (1997).
3. Wolff, B., Sanglier, J.-J., Wang, Y., Leptomycin B is an inhibitor of nuclear export: inhibition of nucleo-cytoplasmic translocation of the human immunodeficiency virus type 1 (HIV-1) Rev protein and Rev-dependent mRNA. *Chem. Biol.* **4**(2), 139-147 (1997).

#### 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, 12/15/2022

#### 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