

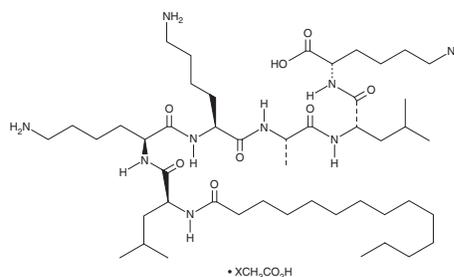
PRODUCT INFORMATION



Myristoyl Hexapeptide-16 (acetate)

Item No. 27549

Formal Name: N-(1-oxotetradecyl)-L-leucyl-L-lysyl-L-lysyl-L-alanyl-L-leucyl-L-lysine, acetate
Synonyms: Myr-Leu-Lys-Lys-Ala-Leu-Lys, Myr-LKKALK
MF: C₄₇H₉₁N₉O₈ • XC₂H₄O₂
FW: 910.3
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

Myristoyl hexapeptide-16 (acetate) is supplied as a solid. A stock solution may be made by dissolving the myristoyl hexapeptide-16 (acetate) in the solvent of choice, which should be purged with an inert gas. Myristoyl hexapeptide-16 (acetate) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of myristoyl hexapeptide-16 (acetate) in these solvents is approximately 30 mg/ml.

Myristoyl hexapeptide-16 (acetate) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, myristoyl hexapeptide-16 (acetate) should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Myristoyl hexapeptide-16 (acetate) has a solubility of approximately 0.25 mg/ml in a 1:3 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Myristoyl hexapeptide-16 is a myristoylated form of the hexapeptide Leu-Lys-Lys-Ala-Leu-Lys. An acid form of myristoyl hexapeptide-16 is active against *E. coli*, *P. aeruginosa*, *S. aureus*, and methicillin-resistant *S. aureus* (MRSA; MICs = ≤100 ppm for all).¹

Reference

1. Owen, D. and Fan, L. Polymeric biosurfactants. *Therapeutic Peptides, Inc.* US 8,586,541 B2 (2013).

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