

PRODUCT INFORMATION



Myristoyl Pentapeptide-17 (acetate)

Item No. 10526

Formal Name: N²-(1-oxotetradecyl)-L-lysyl-L-leucyl-L-alanyl-L-lysyl-L-lysineamide, acetate

Synonyms: Myr-KLAKK-NH₂, Myr-KLAKK-amide, Myr-Lys-Leu-Ala-Lys-Lys-NH₂

MF: C₄₁H₈₁N₉O₆ • XC₂H₄O₂

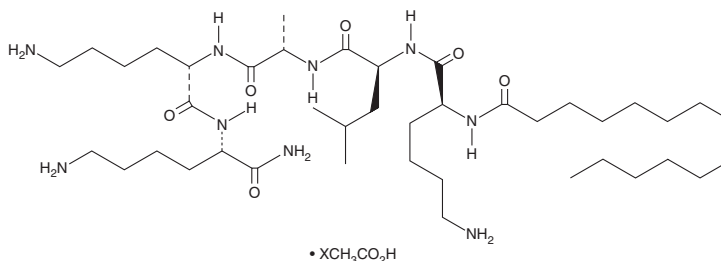
FW: 796.1

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 pentapeptide-17 (acetate) is supplied as a solid. A stock solution may be made by dissolving the myristoyl pentapeptide-17 (acetate) in the solvent of choice, which should be purged with an inert gas. Myristoyl pentapeptide-17 (acetate) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of myristoyl pentapeptide-17 (acetate) in these solvents is approximately 30 mg/ml.

Myristoyl pentapeptide-17 (acetate) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, myristoyl pentapeptide-17 (acetate) should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Myristoyl pentapeptide-17 (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 pentapeptide-17 is a myristoylated form of the pentapeptide Lys-Leu-Ala-Lys-Lys-NH₂. It is active against *E. coli* (MIC = 10 ppm).¹

References

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, 11/08/2022

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