

PRODUCT INFORMATION

Myristoyl Hexapeptide-16 (hydrochloride)

Item No. 40513

Formal Name: tetradecanoyl-L-leucyl-L-lysyl-L-lysyl-L-alanyl-L-leucyl-L-lysine, hydrochloride

Synonym: Myr-Leu-Lys-Lys-Ala-Leu-Lys

Peptide Sequence: Myr-LKKALK

MF: C₄₇H₉₁N₉O₈ • XHCl

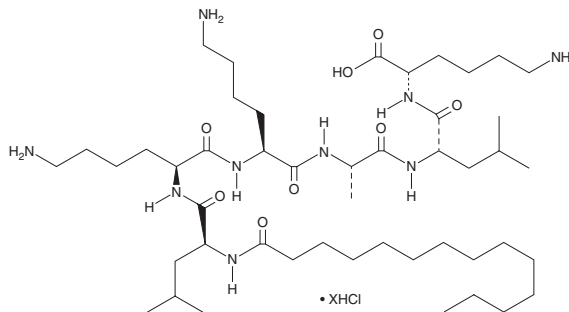
FW: 910.3

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 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the myristoyl hexapeptide-16 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Myristoyl hexapeptide-16 (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and acetonitrile. It is also soluble in water. 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

- 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, 04/10/2024

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