

PRODUCT INFORMATION



LL-37 (human) (scrambled) (trifluoroacetate salt)

Item No. 37503

Peptide Sequence: H-GLKLRFEFSKIKGEFLKTPEVRFDRDIKLDNRISVQR-OH

Synonyms: Cathelicidin, CAP-18, FALL-39, hCAP-18

MF: C₂₀₅H₃₄₀N₆₀O₅₃ • XCF₃COOH

FW: 4,493.3

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.

H—Gly—Leu—Lys—Leu—Arg—Phe—Glu—Phe—Ser—Lys—
Ile—Lys—Gly—Glu—Phe—Leu—Lys—Thr—Pro—Glu—
Val—Arg—Phe—Arg—Asp—Ile—Lys—Leu—Lys—Asp—
Asn—Arg—Ile—Ser—Val—Gln—Arg—OH
• XCF₃COOH

Laboratory Procedures

LL-37 (human) (scrambled) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the LL-37 (human) (scrambled) (trifluoroacetate salt) in water. We do not recommend storing the aqueous solution for more than one day.

Description

LL-37 (human) (scrambled) is a synthetic peptide with identical amino acid composition to LL-37 (human) (Item No. 24461). It has been used as a negative control for LL-37 (human) in various antimicrobial assays.¹

Reference

1. Sheehan, G., Bergsson, G., McElvaney, N.G., *et al.* The human cathelicidin antimicrobial peptide LL-37 promotes the growth of the pulmonary pathogen *Aspergillus fumigatus*. *Infect. Immun.* **86(7)**, e0097-18 (2018).

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, 07/06/2023

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