

PRODUCT INFORMATION



Biotin-PEG₄-LL-37 (human) (trifluoroacetate salt)

Item No. 36996

Peptide Sequence: Biotin-PEG₄-LLGDFFRKSKEKIGKE
FKRIVQRIKDFLRNLVPRTES-OH

MF: C₂₂₆H₃₇₆N₆₄O₅₉S • CF₃COOH

FW: 4,965.9

Purity: ≥90%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

Biotin-PEG₄-Leu-Leu-Gly-Asp-Phe-Phe-Arg-Lys-Ser-Lys-

Glu-Lys-Ile-Gly-Lys-Glu-Phe-Lys-Arg-Ile-

Val-Gln-Arg-Ile-Lys-Asp-Phe-Leu-Arg-Asn-

Leu-Val-Pro-Arg-Thr-Glu-Ser-OH

• XCF₃COOH

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

Laboratory Procedures

Biotin-PEG₄-LL-37 (human) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the biotin-PEG₄-LL-37 (human) (trifluoroacetate salt) in water. We do not recommend storing the aqueous solution for more than one day.

Description

Biotin-PEG₄-LL-37 is a biotinylated and pegylated form of the antimicrobial peptide LL-37 (Item No. 24461).

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, 09/27/2022

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