

PRODUCT INFORMATION



Cys-Pexiganan (trifluoroacetate salt)

Item No. 36987

Peptide Sequence: CGIGKFLKKAKKFGKAFVKILKK-NH₂
MF: C₁₂₅H₂₁₅N₃₃O₂₃S • XCF₃COOH
FW: 2,580.3
Purity: ≥90%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years

H—Cys—Gly—Ile—Gly—Lys—Phe—Leu—Lys—Lys—Ala—
Lys—Lys—Phe—Gly—Lys—Ala—Phe—Val—Lys—Ile—
Leu—Lys—Lys—NH₂
• XCF₃COOH

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

Laboratory Procedures

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

Description

Cys-pexiganan is an antimicrobial peptide and a derivative of pexiganan (Item No. 36924) with an N-terminal cysteine.¹ It is active against *P. aeruginosa* and *S. aureus* (MICs = 64 and 16 µg/ml, respectively). Conjugation of Cys-pexiganan to cellulose acetate-containing polyvinyl alcohol electrospun mats accelerates clotting of isolated recalcified human plasma.

Reference

1. Teixeira, M.A., Antunes, J.C., Seabra, C.L., *et al.* Tiger 17 and pexiganan as antimicrobial and hemostatic boosters of cellulose acetate-containing poly(vinyl alcohol) electrospun mats for potential wound care purposes. *Int. J. Biol. Macromol.* **209(Pt A)**, 1526-1541 (2002).

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, 09/26/2023

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