

PRODUCT INFORMATION



Tiger17 (trifluoroacetate salt)

Item No. 37473

Formal Name: (6S,9R,14R,17S,19aS,25S,27aS,33S,35aS)-N-((S)-1-amino-3-(1H-imidazol-5-yl)-1-oxopropan-2-yl)-9-((S)-2-amino-3-(1H-indol-3-yl)propanamido)-6,25,33-tris(4-aminobutyl)-17-(3-guanidinopropyl)-5,8,16,19,24,27,32,35-octaaxodotriacontahydro-13H-tripyrrolo[2,1-j:2',1'-p:2'',1''-v][1,2]dithia[5,8,11,14,17,20,23,26]octaazacyclononacosine-14-carboxamide, trifluoroacetate salt

Synonym: H-Trp-Cys-Lys-Pro-Lys-Pro-Lys-Pro-Arg-Cys-His-NH₂

Peptide Sequence: WCKPKPKPRCH-NH₂

MF: C₆₂H₉₇N₂₁O₁₁S₂ • XCF₃COOH

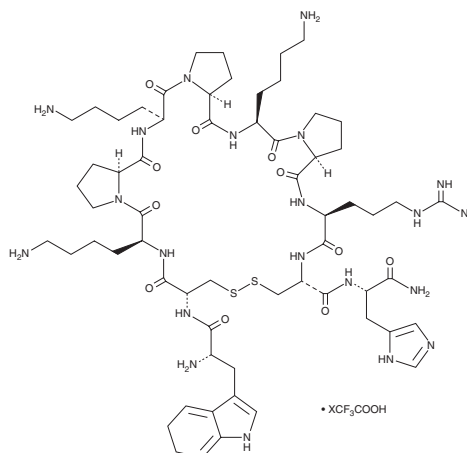
FW: 1,376.7

Purity: ≥98%

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

Tiger17 (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the tiger17 (trifluoroacetate salt) in water. We do not recommend storing the aqueous solution for more than one day.

Description

Tiger17 is a cyclic peptide and derivative of tigerinin-RC1 and tigerinin-RC2, which are antimicrobial peptides found in crab-eating frog (*F. cancrivora*) skin secretions.¹ It increases the proliferation of HaCaT keratinocytes and human skin fibroblasts when used at concentrations ranging from 5 to 20 µg/ml and increases HaCaT cell migration at 20 µg/ml. Tiger17 increases the recruitment of, and the secretion of TGF-β1 and IL-6 from, RAW 264.7 macrophages. Topical application of tiger17 (20 µg/ml) increases the rate of wound closure and re-epithelialization in a mouse model of full-thickness dermal wounds.

Reference

1. Tang, J., Liu, H., Gao, C., *et al.* A small peptide with potential ability to promote wound healing. *PLoS One* **9**(3), e92082 (2014).

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, 10/16/2023

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