

PRODUCT INFORMATION



SAAP 148 (trifluoroacetate salt)

Item No. 37517

Formal Name: (S)-N1-(((S)-1-(((S)-6-amino-1-(((S)-7-amino-1-((S)-1-(((S)-1-(((S)-1-amino-5-guanidino-1-oxopentan-2-yl)amino)-3-methyl-1-oxobutan-2-yl)amino)pyrrolidin-2-yl)-1,2-dioxoheptan-3-yl)amino)-1-oxohexan-2-yl)amino)-4-methyl-1-oxopentan-2-yl)-2-(((2S,5S,8S,11S,14S,17S,20S,23S,26S,29S,32S,35S,38S,41S,44S,47S,50S)-5,38-bis((1H-indol-3-yl)methyl)-14,23,35,47-tetrakis(4-aminobutyl)-26-benzyl-2,11,32,44-tetrakis(3-guanidinopropyl)-8-(4-hydroxybenzyl)-17,20,50-triisobutyl-29,41-diisopropyl-4,7,10,13,16,19,22,25,28,31,34,37,40,43,46,49,52-heptadecaoxo-3,6,9,12,15,18,21,24,27,30,33,36,39,42,45,48,51-heptadecaazatripentacontanamido)pentanediamide, trifluoroacetate salt

Ac—Leu—Lys—Arg—Val—Trp—Lys—Arg—Val—Phe—Lys—
Leu—Leu—Lys—Arg—Tyr—Trp—Arg—Gln—Leu—Lys—
Lys—Pro—Val—Arg—NH₂
• XCF₃COOH

Synonym: Synthetic Antimicrobial and Antibiofilm Peptide 148

Peptide Sequence: Ac-LKRVWKR VFKLLKRYWRQLKKPVR-NH₂

MF: C₁₅₇H₂₆₁N₄₉O₂₇ • XCF₃COOH

FW: 3,267.1

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

SAAP 148 (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the SAAP 148 (trifluoroacetate salt) in water. We do not recommend storing the aqueous solution for more than one day.

Description

SAAP 148 is an antimicrobial peptide.¹ It is active against several multidrug-resistant bacteria *in vitro*. *In vivo*, SAAP 148 (0.125-2% w/w) decreases the number of skin colony-forming units (CFUs) in mouse models of *A. baumannii* or methicillin-resistant *S. aureus* (MRSA) infections.

Reference

1. de Breij, A., Riool, M., Cordfunke, R.A., *et al.* The antimicrobial peptide SAAP-148 combats drug-resistant bacteria and biofilms. *Sci. Transl. Med.* **10(423)**, eaan4044 (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, 10/12/2023

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