

PRODUCT INFORMATION



SAP15 (trifluoroacetate salt)

Item No. 38165

Formal Name: glycyL-L-lysyl-L-cysteinyl-L-seryl-L-threonyl-L-arginylglycyL-L-arginyl-L-lysyl-L-cysteinyl-L-cysteinyl-L-arginyl-L-arginyl-L-lysyl-L-lysine, trifluoroacetate salt

Synonym: Synthetic Anti-inflammatory Peptide 15 H-Gly-Lys-Cys-Ser-Thr-Arg-Gly-Arg-Lys-Cys-Cys-Arg-Arg-Lys-Lys-OH

Peptide Sequence: GKCTRGRKCCRRKK-OH

MF: $C_{68}H_{131}N_{31}O_{18}S_3 \cdot XCF_3COOH$

FW: 1,767.2 • XCF_3COOH

Purity: $\geq 90\%$

Supplied as: A solid

Storage: $-20^\circ C$

Stability: ≥ 4 years

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

Laboratory Procedures

SAP15 (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the SAP15 (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. SAP15 (trifluoroacetate salt) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of SAP15 (trifluoroacetate salt) in these solvents is approximately 0.5, 25, and 10 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of SAP15 (trifluoroacetate salt) can be prepared by directly dissolving the solid in aqueous buffers. The solubility of SAP15 (trifluoroacetate salt) in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

SAP15 is a cell-penetrating peptide consisting of the C-terminal amino acids 52-67 of human β -defensin-3 (Item No. 24578).¹

Reference

1. Lee, J.-Y., Lee, D.W., Jo, B.S., *et al.* Engineered synthetic cell penetrating peptide with intracellular anti-inflammatory bioactivity: An *in vitro* and *in vivo* study. *J. Biomed. Mater. Res. A* **109(10)**, 2001-2016 (2021).

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, 11/20/2023

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