

PRODUCT INFORMATION



SPACE Peptide (trifluoroacetate salt)

Item No. 45602

Formal Name: cyclic (2→10) disulfide, L-alanyl-L-cysteinyl-L-threonylglycyl-L-seryl-L-threonyl-L-glutaminyl-L-histidyl-L-glutaminyl-L-cysteinyl-glycine, trifluoroacetate salt

Synonym: Skin-penetrating and Cell-entering Peptide

Peptide Sequence: Ac-A[CTGSTQHQC]G-OH

MF: C₄₀H₆₃N₁₅O₁₇S₂ • XCF₃COOH

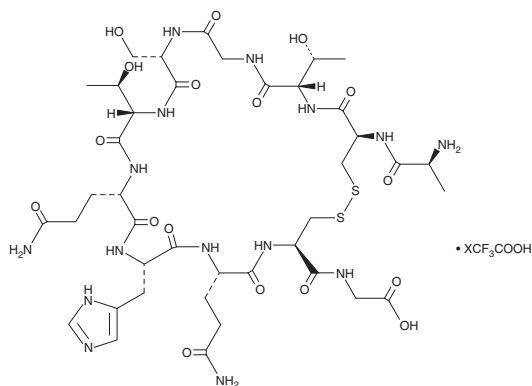
FW: 1,090.2

Purity: ≥95%

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

SPACE peptide (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the SPACE peptide (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. SPACE peptide (trifluoroacetate salt) is slightly soluble (0.1-1 mg/ml) in DMSO.

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 SPACE peptide (trifluoroacetate salt) can be prepared by directly dissolving the solid in aqueous buffers. SPACE peptide (trifluoroacetate salt) is sparingly soluble (1-10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

SPACE peptide is a skin-penetrating and cell-entering (SPACE) peptide.¹ It is internalized in keratinocytes, fibroblasts, human umbilical vein endothelial cells (HUVECs), and MDA-MB-231 breast cancer cells via macropinocytosis, a form of endocytosis. SPACE peptide-containing conjugates have been used to deliver proteins and siRNA *in vitro* and *in vivo*.

Reference

1. Hsu, T. and Mitragotri, S. Delivery of siRNA and other macromolecules into skin and cells using a peptide enhancer. *Proc. Natl. Acad. Sci. U S A* **108**(38), 15816-15821 (2011).

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, 06/09/2026

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