

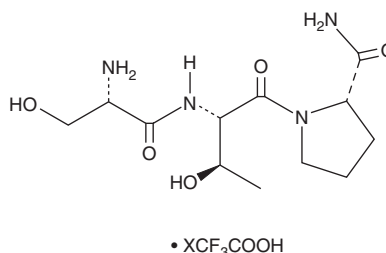
PRODUCT INFORMATION



Tripeptide-32 (trifluoroacetate salt)

Item No. 45955

Formal Name: L-seryl-L-threonyl-L-prolinamide, trifluoroacetate salt
Synonym: Ser-Thr-Pro-NH₂
Peptide Sequence: STP-NH₂
MF: C₁₂H₂₂N₄O₅ • XCF₃COOH
FW: 302.3
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

Tripeptide-32 (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the tripeptide-32 (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Tripeptide-32 (trifluoroacetate salt) is sparingly soluble (1-10 mg/ml) in ethanol and 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 tripeptide-32 (trifluoroacetate salt) can be prepared by directly dissolving the solid in aqueous buffers. Tripeptide-32 (trifluoroacetate salt) is soluble (≥10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

Tripeptide-32 is a cosmeceutical peptide.¹ Formulations containing tripeptide-32 have been used in anti-aging cosmetics.

Reference

1. He, Q., Liao, Y., Wu, Y., et al. Bioactive oligopeptides and the application in skin regeneration and rejuvenation. *J. Appl. Biomater. Funct. Mater.* (2025).

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, 07/20/2026

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