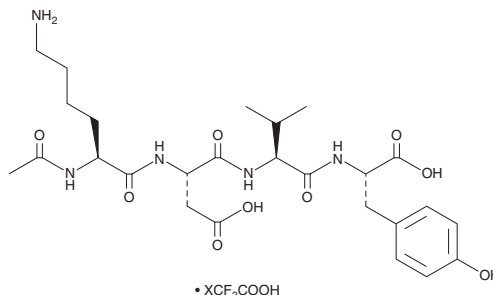


PRODUCT INFORMATION

Acetyl Tetrapeptide-2 (trifluoroacetate salt)

Item No. 45935

Formal Name: N²-acetyl-L-lysyl-L-α-aspartyl-L-valyl-L-tyrosine, trifluoroacetate salt
Synonym: Acetyl-L-Lys-L-Asp-L-Val-L-Tyr-OH
Peptide Sequence: Ac-KDVY-OH
MF: C₂₆H₃₉N₅O₉ • XCF₃COOH
FW: 565.6
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

Acetyl tetrapeptide-2 (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the acetyl tetrapeptide-2 (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Acetyl tetrapeptide-2 (trifluoroacetate salt) is soluble (≥10 mg/ml) in ethanol and sparingly soluble (1-10 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 acetyl tetrapeptide-2 (trifluoroacetate salt) can be prepared by directly dissolving the solid in aqueous buffers. Acetyl tetrapeptide-2 (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

Acetyl tetrapeptide-2 is a peptide derivative of thymopoietin II that corresponds to amino acids 33-36 of the human sequence.¹ It increases the stiffness of HaCaT keratinocytes at 0.5 µg/ml.² Formulations containing acetyl tetrapeptide-2 have been used in cosmetics as skin firming agents.

References

1. Harris, C.A., Andryuk, P.J., Cline, S., *et al.* Three distinct human thymopoietins are derived from alternatively spliced mRNAs. *Proc. Natl. Acad. Sci. U.S.A.* **91**(14), 6283-6287 (1994).
2. Kobiela, T., Milner-Krawczyk, M., Pasikowska-Piwko, M., *et al.* The effect of anti-aging peptides on mechanical and biological properties of HaCaT keratinocytes. *Int. J. Pept. Res. Ther.* **24**(4), 577-587 (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, 06/30/2026

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