

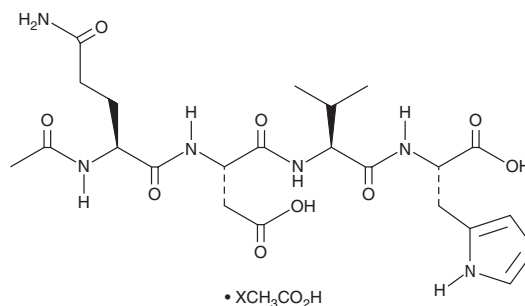
# PRODUCT INFORMATION



## Acetyl Tetrapeptide-9 (acetate)

Item No. 35181

**Formal Name:** N<sup>2</sup>-acetyl-L-glutamyl-L-α-aspartyl-L-valyl-L-histidine, acetate  
**Synonyms:** AcTP1, N-acetyl-Gln-Asp-Val-His  
**MF:** C<sub>22</sub>H<sub>33</sub>N<sub>7</sub>O<sub>9</sub> • XC<sub>2</sub>H<sub>4</sub>O<sub>2</sub>  
**FW:** 539.5  
**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

Acetyl tetrapeptide-9 (acetate) is supplied as a solid. Aqueous solutions of acetyl tetrapeptide-9 (acetate) can be prepared by directly dissolving the acetyl tetrapeptide-9 (acetate) in aqueous buffers. The solubility of acetyl tetrapeptide-9 (acetate) in PBS (pH 7.2) is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Acetyl tetrapeptide-9 is a synthetic signal peptide.<sup>1</sup> It stimulates the synthesis of lumican *in vitro*.<sup>2</sup>

### References

1. Gorouhi, F. and Maibach, H.I. Role of topical peptides in preventing or treating aged skin. *Int. J. Cosmet. Sci.* **31(5)**, 327-345 (2009).
2. Pauly, G., Contet-Audonneau, J.-L., Moussou, P., et al. Small proteoglycans in the skin: New targets in the fight against skin aging. *Int. J. Cosmet. Sci.* **31(2)**, 154 (2009).

#### 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, 11/21/2022

#### 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