

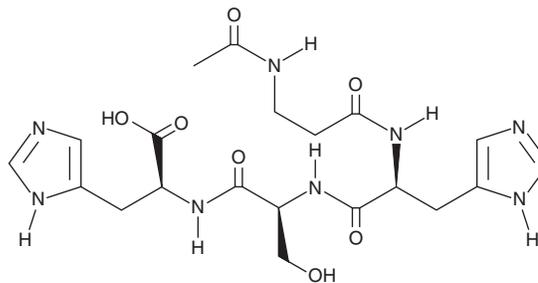
# PRODUCT INFORMATION



## Acetyl Tetrapeptide-5

Item No. 27152

**CAS Registry No.:** 820959-17-9  
**Formal Name:** N-acetyl-β-alanyl-L-histidyl-L-seryl-L-histidine  
**MF:** C<sub>20</sub>H<sub>28</sub>N<sub>8</sub>O<sub>7</sub>  
**FW:** 492.5  
**Purity:** ≥98%  
**Supplied as:** A crystalline 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-5 is supplied as a crystalline solid. Aqueous solutions of acetyl tetrapeptide-5 can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of acetyl tetrapeptide-5 in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Acetyl tetrapeptide-5 is a cosmeceutical matricin peptide.<sup>1</sup>

### Reference

- Schagen, S.K. Topical peptide treatments with effective anti-aging results. *Cosmetics* **4**(2), (2017).

#### 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, 12/13/2022

#### 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