

# PRODUCT INFORMATION



## Palmitoyl Tripeptide-38 (acetate)

Item No. 27564

**Formal Name:** ((S)-2-((S)-6-amino-2-palmitamido-hexanamido)-4-(methylsulfonyl)butanoyl)-L-lysine, acetic acid

**Synonyms:** Pal-KMO<sub>2</sub>K-OH, Pal-Lys-Met(O<sub>2</sub>)-Lys-OH

**MF:** C<sub>33</sub>H<sub>65</sub>N<sub>5</sub>O<sub>7</sub>S • XC<sub>2</sub>H<sub>4</sub>O<sub>2</sub>

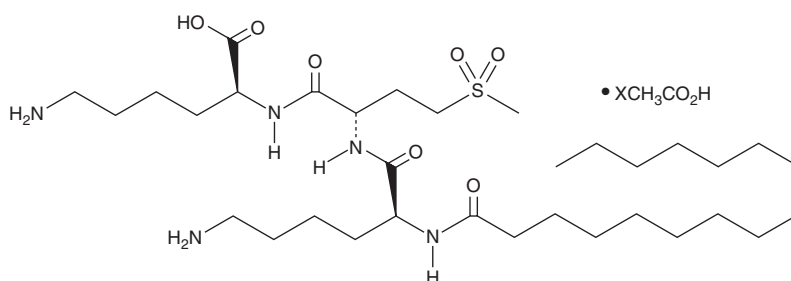
**FW:** 676.0

**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

Palmitoyl tripeptide-38 (acetate) is supplied as a solid. A stock solution may be made by dissolving the palmitoyl tripeptide-38 (acetate) in the solvent of choice, which should be purged with an inert gas. Palmitoyl tripeptide-38 (acetate) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of palmitoyl tripeptide-38 (acetate) in these solvents is approximately 30 mg/ml.

Palmitoyl tripeptide-38 (acetate) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, palmitoyl tripeptide-38 (acetate) should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Palmitoyl tripeptide-38 (acetate) has a solubility of approximately 0.25 mg/ml in a 1:3 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Palmitoyl tripeptide-38 is a palmitoylated form of the tripeptide Lys-Met(O<sub>2</sub>)-Lys-OH.<sup>1</sup> It increases the levels of collagen I, collagen II, fibronectin, and hyaluronic acid by 187, 97, 109, and 55%, respectively, produced by isolated human dermal fibroblasts when used at a concentration of 6 ppm.<sup>2</sup>

### References

- Schagen, S.K. Topical peptide treatments with effective anti-aging results. *Cosmetics* **4**(2), (2017).
- Fournial, A., Mondon, P., and Peschard, O. KXX-Peptide compositions for cosmetic and dermo-pharmaceutic uses. *Sederma Le Perray en Yvelines*. EP2382230A2 (2010).

#### 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, 12/19/2022

#### 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