

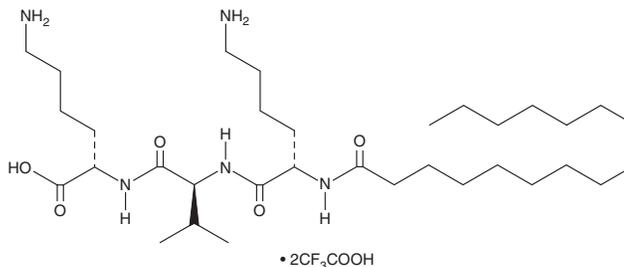
# PRODUCT INFORMATION



## Palmitoyl Tripeptide-5 (trifluoroacetate salt)

Item No. 27565

**CAS Registry No.:** 623172-56-5  
**Formal Name:** N<sup>2</sup>-(1-oxohexadecyl)-L-lysyl-L-valyl-L-lysine, 2,2,2-trifluoroacetate (1:2)  
**MF:** C<sub>33</sub>H<sub>65</sub>N<sub>5</sub>O<sub>5</sub> • 2CF<sub>3</sub>COOH  
**FW:** 840.0  
**Purity:** ≥95%  
**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

Palmitoyl tripeptide-5 (trifluoroacetate salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the palmitoyl tripeptide-5 (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Palmitoyl tripeptide-5 (trifluoroacetate salt) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of palmitoyl tripeptide-5 (trifluoroacetate salt) in these solvents is approximately 30 mg/ml.

Palmitoyl tripeptide-5 (trifluoroacetate salt) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, palmitoyl tripeptide-5 (trifluoroacetate salt) should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Palmitoyl tripeptide-5 (trifluoroacetate salt) 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-5 is a cosmeceutical peptide.<sup>1</sup>

### Reference

1. Bidamant, F., Imflid, D., Joller, P., *et al.* Novel compositions and their use. *DSM IP Assets B.V. WO 2009/124971 A2* (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, 12/07/2022

#### 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