

# PRODUCT INFORMATION



## Hexapeptide-11 (acetate)

Item No. 27155

**Formal Name:** L-phenylalanyl-L-valyl-L-alanyl-L-prolyl-L-phenylalanyl-L-proline, acetic acid

**Synonyms:** FVAPFP, Phe-Val-Ala-Pro-Phe-Pro

**MF:** C<sub>36</sub>H<sub>48</sub>N<sub>6</sub>O<sub>7</sub> • XC<sub>2</sub>H<sub>4</sub>O<sub>2</sub>

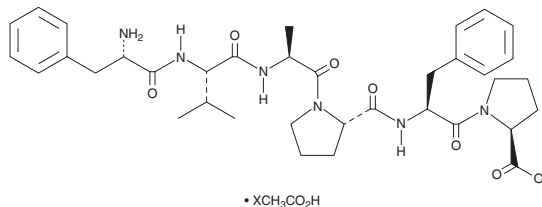
**FW:** 676.8

**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

Hexapeptide-11 (acetate) is supplied as a crystalline solid. A stock solution may be made by dissolving the hexapeptide-11 (acetate) in the solvent of choice, which should be purged with an inert gas. Hexapeptide-11 (acetate) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of hexapeptide-11 (acetate) in these solvents is approximately 30 mg/ml.

### Description

Hexapeptide-11 is a synthetic hexapeptide.<sup>1</sup> Hexapeptide-11 (5% v/v) increases the expression of the antioxidant genes *Nrf2*, *Keap1*, *NQO1*, and *TXNRD1* and decreases hydrogen peroxide-induced senescence of IMR-90 human lung fibroblasts *in vitro*. Hexapeptide-11 (5% v/v) increases the activity of matrix metalloproteinase-2 (MMP-2) and MMP-9 and the proteasome peptidases LLVY/β5 and LLE/β1 in IMR-90 cells *in vitro* and reduces the migration of IMR-90 cells in a scratch wound assay.

### Reference

1. Sklirou, A.D., Ralli, M., Dominguez, M., *et al.* Hexapeptide-11 is a novel modulator of the proteostasis network in human diploid fibroblasts. *Redox Biol.* **5**, 205-215 (2015).

#### 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, 10/20/2022

#### 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