

PRODUCT INFORMATION



Acetyl Hexapeptide-49 (trifluoroacetate salt)

Item No. 45626

Formal Name: (S)-2-((2S,5S,8S,11S,14S)-2-((1H-imidazol-4-yl)methyl)-8-((1H-indol-3-yl)methyl)-5,11,14-tribenzyl-4,7,10,13,16-pentaoxo-3,6,9,12,15-pentaazaheptadecanamido)-3-methylbutanamide, trifluoroacetate salt

Synonym: Ac-Phe-Phe-Trp-Phe-His-Val-NH₂

Peptide Sequence: Ac-FFWFHV-NH₂

MF: C₅₁H₅₈N₁₀O₇ • XCF₃COOH

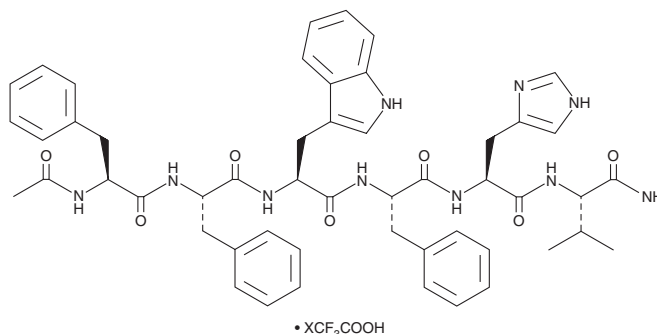
FW: 923.1

Purity: ≥98%

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 hexapeptide-49 (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the acetyl hexapeptide-49 (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Acetyl hexapeptide-49 (trifluoroacetate salt) is sparingly soluble (1-10 mg/ml) in acetonitrile:water (1:1).

Description

Acetyl hexapeptide-49 is an antagonist of proteinase-activated receptor 2 (PAR2).^{1,2} It inhibits PAR2 agonist-induced calcium release and calcitonin gene-related peptide (CGRP) and IL-6 production in human keratinocytes.¹ Acetyl hexapeptide-49 also induces proliferation in human keratinocytes and has photoprotective activity in irradiated human dermal fibroblasts. Formulations containing acetyl hexapeptide-49 have been used in cosmetics.

References

1. Yau, M.-K., Lim, J., Liu, L., *et al.* Protease activated receptor 2 (PAR2) modulators: A patent review (2010-2015). *Expert Opin. Ther. Pat.* **26**(4), 471-483 (2016).
2. Resende, D.I.S.P., Ferreira, M.S., Sousa-Lobo, J.M., *et al.* Usage of synthetic peptides in cosmetics for sensitive skin. *Pharmaceuticals (Basel)* **14**(8), 702 (2021).

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, 06/02/2026

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