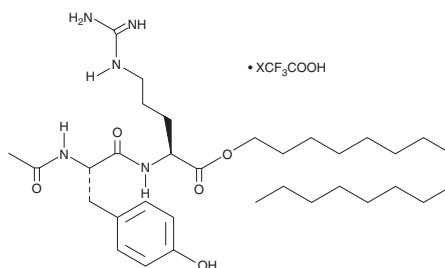


# PRODUCT INFORMATION

## Acetyl Dipeptide 1 cetyl ester (trifluoroacetate salt)

Item No. 34949

**Formal Name:** N-acetyl-L-tyrosyl-L-arginine, hexadecyl ester, trifluoroacetate salt  
**MF:**  $C_{33}H_{57}N_5O_5 \cdot XCF_3COOH$   
**FW:** 603.8  
**Purity:**  $\geq 98\%$   
**UV/Vis.:**  $\lambda_{max}$ : 225 nm  
**Supplied as:** A solid  
**Storage:**  $-20^\circ C$   
**Stability:**  $\geq 4$  years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

Acetyl dipeptide 1 cetyl ester (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the acetyl dipeptide 1 cetyl ester (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Acetyl dipeptide 1 cetyl ester (trifluoroacetate salt) is soluble in DMSO.

### Description

Acetyl dipeptide 1 cetyl ester is a cosmeceutical peptide.<sup>1</sup> It increases expression of the genes encoding the skin barrier function proteins AQP3, FLG, caspase-14, and cytokeratin 10 in primary human keratinocytes. Acetyl dipeptide 1 cetyl ester reduces glyoxal-induced formation of advanced glycation end products (AGEs) in primary human fibroblasts.

### Reference

1. Khmaladze, I., Österlund, C., Smiljanic, S., *et al.* A novel multifunctional skin care formulation with a unique blend of antipollution, brightening and antiaging active complexes. *J. Cosmet. Dermatol.* **19**(6), 1415-1425 (2020)

#### 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, 10/12/2023

#### 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