

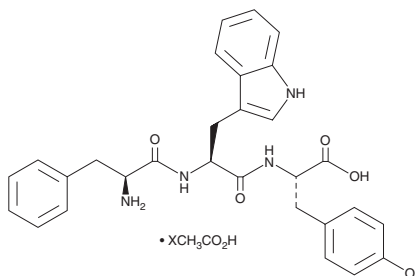
# PRODUCT INFORMATION



## Tripeptide-41 (acetate)

Item No. 45598

**Formal Name:** L-phenylalanyl-L-tryptophyl-L-tyrosine, acetate  
**Synonym:** H-Phe-Trp-Tyr-OH  
**Peptide Sequence:** H-FWY-OH  
**MF:**  $C_{29}H_{30}N_4O_5 \cdot XC_2H_4O_2$   
**FW:** 514.6  
**Purity:**  $\geq 98\%$   
**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

Tripeptide-41 (acetate) is supplied as a solid. A stock solution may be made by dissolving the tripeptide-41 (acetate) in the solvent of choice, which should be purged with an inert gas. Tripeptide-41 (acetate) is soluble ( $\geq 10$  mg/ml) in ethanol. Tripeptide-41 (acetate) is sparingly soluble (1-10 mg/ml) in DMSO.

### Description

Tripeptide-41 is a cosmeceutical peptide.<sup>1</sup>

### Reference

1. Marini, J.L. and Saxena, S.J. Anti-cellulite cream. *Jan Marini Skin Research* (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, 06/03/2026

#### 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](http://WWW.CAYMANCHEM.COM)