

PRODUCT INFORMATION



Decapeptide-12 (acetate)

Item No. 35400

Formal Name: L-tyrosyl-L-arginyl-L-seryl-L-arginyl-L-lysyl-L-tyrosyl-L-seryl-L-seryl-L-tryptophyl-L-tyrosine, acetate

Synonym: YRSRKYSSWY

MF: C₆₅H₉₀N₁₈O₁₇ • XC₂H₄O₂

FW: 1,395.5

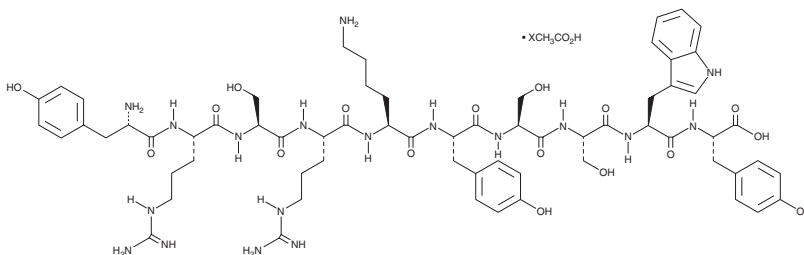
Purity: ≥95%

UV/Vis.: λ_{max}: 221 nm

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

Decapeptide-12 (acetate) is supplied as a solid. A stock solution may be made by dissolving the decapeptide-12 (acetate) in water. We do not recommend storing the aqueous solution for more than one day.

Description

Decapeptide-12 is a tyrosinase inhibitor and cosmeceutical peptide.¹ It inhibits mushroom tyrosinase (IC₅₀ = 40 μM), as well as human tyrosinase by 35% when used at a concentration of 100 μM.

Reference

1. Hariri, R.J., Saeedi, M., and Akbarzadeh, T. Naturally occurring and synthetic peptides: Efficient tyrosinase inhibitors. *J. Pept. Sci.* **27(7)**, e3329 (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, 12/12/2022

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