

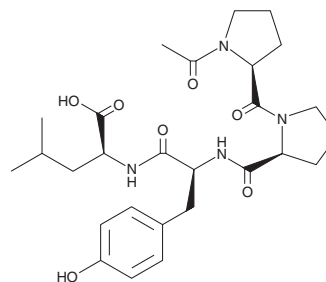
PRODUCT INFORMATION



Acetyl Tetrapeptide-11

Item No. 42985

CAS Registry No.: 928006-88-6
Formal Name: 1-acetyl-L-prolyl-L-prolyl-L-tyrosyl-L-leucine
Synonym: Ac-Pro-Pro-Tyr-Leu-OH
Peptide Sequence: Ac-PPYL-OH
MF: C₂₇H₃₈N₄O₇
FW: 530.6
Purity: ≥95%
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 tetrapeptide-11 is supplied as a solid. A stock solution may be made by dissolving the acetyl tetrapeptide-11 in the solvent of choice, which should be purged with an inert gas. Acetyl tetrapeptide-11 is soluble (≥10 mg/ml) in organic solvents such as ethanol and DMSO.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of acetyl tetrapeptide-11 can be prepared by directly dissolving the solid in aqueous buffers. Acetyl tetrapeptide-11 is sparingly soluble (1-10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

Acetyl tetrapeptide-11 is a peptide inducer of syndecan-1 (SYND1) and collagen XVII.¹ It increases SYND1 and collagen XVII levels in primary human keratinocytes when used at a concentration of 2.6 µg/ml. Formulations containing acetyl tetrapeptide-11 have been used in skin care cosmetics.

Reference

1. Rodrigues, A.-L., Benoit, I., Clarius, T., *et al.* Counteracting ageing phenomena by new pure tetrapeptides with targeted efficacy. *J. Appl. Cosmetol.* **27(2)**, 63-71 (2008).

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, 07/07/2025

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