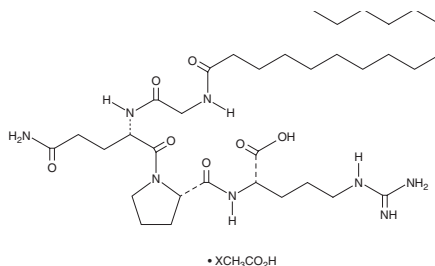


PRODUCT INFORMATION

Palmitoyl Tetrapeptide-7 (acetate)

Item No. 27550

Formal Name: N-(1-oxohexadecyl)glycyl-L-glutamyl-L-prolyl-L-arginine
Synonyms: N-Palmitoylargin, Pal-Gly-Gln-Pro-Arg-OH
MF: $C_{34}H_{62}N_8O_7 \cdot XC_2H_4O_2$
FW: 694.9
Purity: $\geq 98\%$
Supplied as: A solid
Storage: $-20^\circ C$
Stability: ≥ 4 years



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

Laboratory Procedures

Palmitoyl tetrapeptide-7 (acetate) is supplied as a solid. A stock solution may be made by dissolving the palmitoyl tetrapeptide-7 (acetate) in the solvent of choice, which should be purged with an inert gas. Palmitoyl tetrapeptide-7 (acetate) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of palmitoyl tetrapeptide-7 (acetate) in these solvents is approximately 5, 15, and 20 mg/ml, respectively.

Description

Palmitoyl tetrapeptide-7 is a cosmeceutical peptide that stimulates type I collagen, fibronectin, and hyaluronic acid production in skin.¹ Formulations containing palmitoyl tetrapeptide-7 have been used in the treatment of photoaging, dark circles, and undereye puffiness.

Reference

1. Bucay, V.W. and Day, D. Adjunctive skin care of the brow and periorbital region. *Clin. Plast. Surg.* **40(1)**, 225-236 (2013).

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, 10/10/2022

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