

PRODUCT INFORMATION



Pal-VGVAPG (acetate)

Item No. 35185

Formal Name: N-(1-oxohexadecyl)-L-valylglycyl-L-valyl-L-alanyl-L-prolyl-glycine, acetate

Synonyms: Palmitoyl Hexapeptide-12, Palmitoyl-VGVAPG, Pal-Val-Gly-Val-Ala-Pro-Gly

MF: $C_{38}H_{68}N_6O_8 \cdot XC_2H_4O_2$

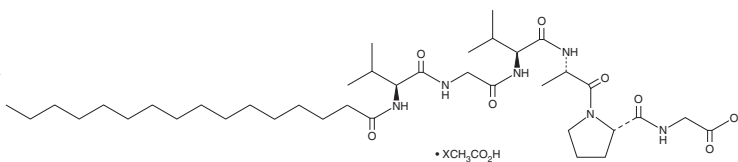
FW: 737.0

Purity: $\geq 98\%$

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

Pal-VGVAPG (acetate) is supplied as a solid. A stock solution may be made by dissolving the Pal-VGVAPG (acetate) in the solvent of choice, which should be purged with an inert gas. Pal-VGVAPG (acetate) is soluble in the organic solvent dimethyl formamide at a concentration of approximately 1 mg/ml.

Description

Pal-VGVAPG is a lipidated hexapeptide composed of a fragment of elastin conjugated to palmitic acid (Item No. 10006627).¹ Formulations containing Pal-VGVAPG have been used in cosmeceuticals.

Reference

1. Resende, D.I.S.P., Ferreira, M.S., Sousa-Lobo, J.M., *et al.* Usage of synthetic peptides in cosmetics for sensitive skin. *Pharmaceuticals (Basel)* **14**(8), 702 (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