

PRODUCT INFORMATION



Palmitoyl Tetrapeptide-10 (hydrochloride)

Item No. 45965

Formal Name: N-(1-oxohexadecyl)-L-lysyl-L-threonyl-L-phenylalanyl-L-lysine, hydrochloride

Synonyms: Pal-Lys-Thr-Phe-Lys-OH, Palmitoyl-KTFK, Palmitoyl-Lys-Thr-Phe-Lys

Peptide Sequence: Pal-KTFK-OH

MF: $C_{41}H_{72}N_6O_7 \cdot XHCl$

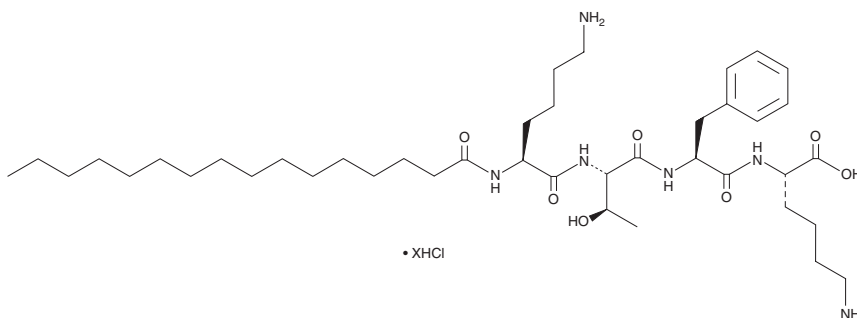
FW: 761.1

Purity: $\geq 95\%$

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-10 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the palmitoyl tetrapeptide-10 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Palmitoyl tetrapeptide-10 (hydrochloride) is soluble (≥ 10 mg/ml) in 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 palmitoyl tetrapeptide-10 (hydrochloride) can be prepared by directly dissolving the solid in aqueous buffers. Palmitoyl tetrapeptide-10 (hydrochloride) is soluble (≥ 10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

Palmitoyl tetrapeptide-10 is a lipidated tetrapeptide composed of a four-amino acid peptide fragment conjugated to palmitic acid (Item Nos. 10006627 | 43561). Formulations containing palmitoyl tetrapeptide-10 have been used in cosmetics.

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, 07/17/2026

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