

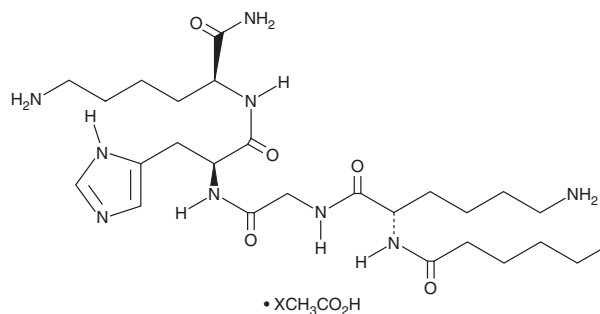
PRODUCT INFORMATION



Caprooyl Tetrapeptide-3 (acetate)

Item No. 27154

CAS Registry No.: 1012317-71-3
Formal Name: N²-(1-oxohexyl)-L-lysylglycyl-L-histidyl-L-lysineamide, acetate
Synonyms: KGHK-Caproic Acid, Lys-Gly-His-Lys-Caproic Acid
MF: C₂₆H₄₇N₉O₅ • XC₂H₄O₂
FW: 565.7
Purity: ≥98%
Supplied as: A crystalline 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

Caprooyl tetrapeptide-3 (acetate) is supplied as a crystalline solid. A stock solution may be made by dissolving the caprooyl tetrapeptide-3 (acetate) in the solvent of choice, which should be purged with an inert gas. Caprooyl tetrapeptide-3 (acetate) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of caprooyl tetrapeptide-3 (acetate) is approximately 30 mg/ml in ethanol and approximately 25 mg/ml in DMSO and DMF.

Caprooyl tetrapeptide-3 (acetate) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, caprooyl tetrapeptide-3 (acetate) should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Caprooyl tetrapeptide-3 (acetate) has a solubility of approximately 0.25 mg/ml in a 1:3 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Caprooyl tetrapeptide-3 is a derivative of tetrapeptide-3 that is linked to caproic acid.¹ It has migration-imparting effects in human adipose-derived stem cells when applied at concentrations of 0.1 and 1 mg/ml.

Reference

1. Tsuda, T. and Yumoto, M. Cosmetic method and dermatological topical agent for use therein, migration imparting agent, and method for screening components for use in cosmetic method for improving skin state. *Rohto Pharmaceutical Co., Ltd. US, US20180161262A1* (2018).

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, 11/07/2022

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