

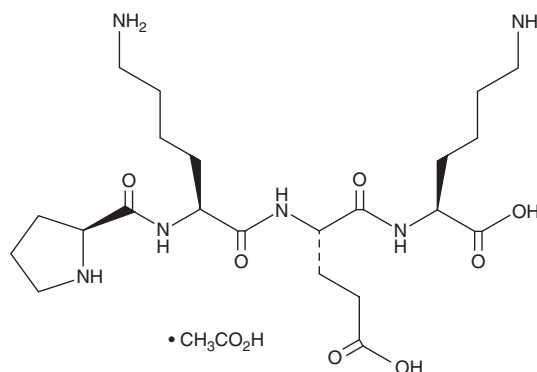
PRODUCT INFORMATION



Tetrapeptide PKEK (acetate)

Item No. 45774

CAS Registry No.: 2763584-89-8
Formal Name: L-prolyl-L-lysyl-L-glutamyl-L-lysine, monoacetate
Synonyms: Pro-Lys-Glu-Lys-OH, Tetrapeptide-30
Peptide Sequence: PKEK-OH
MF: $C_{22}H_{40}N_6O_7 \cdot C_2H_4O_2$
FW: 620.7
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

Tetrapeptide PKEK (acetate) is supplied as a solid. Aqueous solutions of tetrapeptide PKEK (acetate) can be prepared by directly dissolving the solid in aqueous buffers. Tetrapeptide PKEK (acetate) 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

Tetrapeptide PKEK is a bioactive peptide.¹ It reduces UVB-induced expression of proopiomelanocortin (POMC), IL-6, and TNF- α in primary human epidermal keratinocytes at 1 and 10 $\mu\text{g/ml}$ and IL-8 at 10 $\mu\text{g/ml}$.

Reference

1. Marini, A., Farwick, M., Grether-Beck, S., *et al.* Modulation of skin pigmentation by the tetrapeptide PKEK: *In vitro* and *in vivo* evidence for skin whitening effects. *Exp. Dermatol.* **21(2)**, 140-146 (2012).

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, 07/28/2026

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