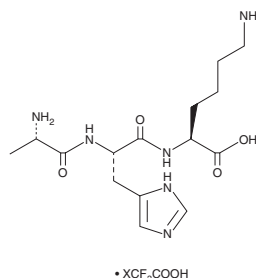


PRODUCT INFORMATION

AHK (trifluoroacetate salt)

Item No. 45938

Formal Name: L-alanyl-L-histidyl-L-lysine, trifluoroacetate salt
Synonym: Ala-His-Lys-OH
Peptide Sequence: AHK-OH
MF: $C_{15}H_{26}N_6O_4 \cdot XCF_3COOH$
FW: 354.4
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

AHK (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the AHK (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. AHK (trifluoroacetate salt) is slightly soluble (0.1-1 mg/ml) in DMSO and acetonitrile:water (1:1).

Description

AHK is a tripeptide that binds to copper to form AHK-Cu, a complex that stimulates hair growth.¹ AHK increases the proliferation of primary human hair follicle dermal papilla cells.²

References

1. Pyo, H.K., Yoo, H.G., Won, C.H., et al. The effect of tripeptide-copper complex on human hair growth *in vitro*. *Arch. Pharm. Res.* **30(7)**, 834-839 (2007).
2. Jeong, J., Park, K.-Y., and Kim, J. Synthesis and biological evaluation of dimeric peptide derivatives as proliferation-stimulating agents in human follicle dermal papilla cells. *Bull. Korean Chem. Soc.* **42(11)**, 1435-1439 (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, 07/15/2026

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