

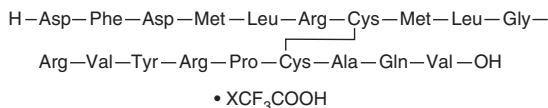
PRODUCT INFORMATION



[Ala¹⁷]-MCH (trifluoroacetate salt)

Item No. 35262

Formal Name: 17-L-alanine-melanin-concentrating hormone (human), trifluoroacetate salt
Synonym: [Ala¹⁷]-Melanin-Concentrating Hormone
MF: C₉₇H₁₅₅N₂₉O₂₆S₄ • XCF₃COOH
FW: 2,271.7
Purity: ≥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

[Ala¹⁷]-MCH (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the [Ala¹⁷]-MCH (trifluoroacetate salt) in water. The solubility of [Ala¹⁷]-MCH (trifluoroacetate salt) in water is approximately 2 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

[Ala¹⁷]-MCH is a synthetic peptide derived from the hypothalamic neuropeptide hormone melanin-concentrating hormone (MCH) and an MCH receptor agonist.¹ It binds to MCH₁ and MCH₂ receptors in CHO cells expressing the human receptors (IC₅₀s = 0.15 and 3.5 nM, respectively) and induces calcium mobilization in HEK293 cells expressing the human receptors (EC₅₀s = 17 and 54 nM, respectively).

Reference

1. Bednarek, M.A., Feighner, S.D., Hreniuk, D.L., *et al.* Short segment of human melanin-concentrating hormone that is sufficient for full activation of human melanin-concentrating hormone receptors 1 and 2. *Biochemistry* **40**(31), 9379-9386 (2001).

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, 11/17/2022

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