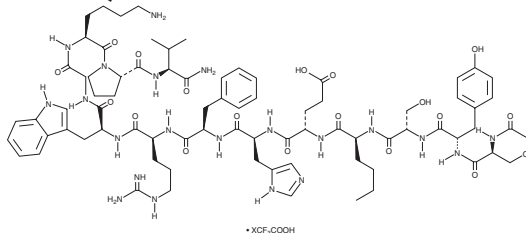


PRODUCT INFORMATION

[Nle⁴,D-Phe⁷]-α-MSH (trifluoroacetate salt)

Item No. 24475

MF: C₇₈H₁₁₁N₂₁O₁₉ • XCF₃COOH
FW: 1,646.9
Purity: ≥95%
Supplied as: A lyophilized powder
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

[Nle⁴,D-Phe⁷]-α-MSH (trifluoroacetate salt) is supplied as a lyophilized powder. A stock solution may be made by dissolving the [Nle⁴,D-Phe⁷]-α-MSH (trifluoroacetate salt) in water. The solubility of [Nle⁴,D-Phe⁷]-α-MSH (trifluoroacetate salt) in water is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

[Nle⁴,D-Phe⁷]-α-MSH is a synthetic analog of α-melanocyte stimulating hormone (α-MSH).¹ It is 26-fold more potent than α-MSH at inducing adenylate cyclase activity in mouse melanoma cells *in vitro*. [Nle⁴,D-Phe⁷]-α-MSH also induces tyrosinase activity in a time-dependent manner in mouse melanoma cells. *In vivo*, [Nle⁴,D-Phe⁷]-α-MSH (1.8 μg/g) induces long-term skin darkening in *R. pipiens* (frogs).² Formulations containing [Nle⁴,D-Phe⁷]-α-MSH have been used to induce melanogenesis in patients with erythropoietic protoporphyria.

References

1. Sawyer, T.K., Sanfilippo, P.J., Hruby, V.J., *et al.* 4-Norleucine, 7-D-phenylalanine-α-melanocyte-stimulating hormone: A highly potent α-melanotropin with ultralong biological activity. *Proc. Nat. Acad. Sci. USA* **77**(10), 5754-5758 (1980).
2. Hadley, M.E., Anderson, B., Heward, C.B., *et al.* Calcium-dependent prolonged effects on melanophores of [4-norleucine, 7-D-phenylalanine]-α-melanotropin. *Science* **213**(4511), 1025-1027 (1981).

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, 12/19/2022

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