

PRODUCT INFORMATION

α-MSH Acid (human, mouse, rat, porcine, bovine, ovine) (trifluoroacetate salt)

Item No. 41463

Formal Name: 13-L-valine- α -melanotropin, trifluoroacetate salt

Synonyms: N-acetyl ACTH (1-13), N-acetyl Adrenocorticotrophic Hormone (1-13), α -Melanocyte-stimulating Hormone Acid

Peptide Sequence: Ac-SYSMEHFRWGKPV-OH

MF: $C_{77}H_{108}N_{20}O_{20}S \bullet XCF_3COOH$

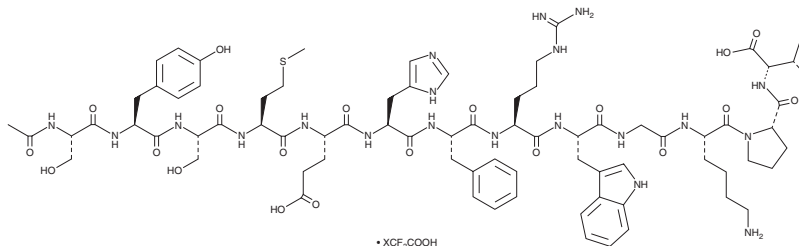
FW: 1,665.9

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

α -MSH acid (human, mouse, rat, porcine, bovine, ovine) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the α -MSH acid (human, mouse, rat, porcine, bovine, ovine) (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. α -MSH acid (human, mouse, rat, porcine, bovine, ovine) (trifluoroacetate salt) is soluble (≥ 10 mg/ml) in the organic solvent DMSO.

Description

α-Melanocyte-stimulating hormone (α-MSH) acid is an acid form of α-MSH (Item No. 29923).¹ It selectively binds to melanocortin receptor 1 (MC1R) over MC3R, MC4R, and MC5R in COS-1 cells expressing the human receptors (K_s = 0.0012, 0.0829, 3.05, and 18.2 μM, respectively).

Reference

1. Schiöth, H.B., Muceniece, R., Larsson, M., *et al.* The melanocortin 1, 3, 4 or 5 receptors do not have a binding epitope for ACTH beyond the sequence of α -MSH. *J. Endocrinol.* 155(1), 73-78 (1997).

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, 10/09/2024

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