

PRODUCT INFORMATION



HS-014 (trifluoroacetate salt)

Item No. 41556

Formal Name: cyclic (1→8)-disulfide N-acetyl-L-cysteinyl-L-α-glutamyl-L-histidyl-3-(2-naphthalenyl)-D-alanyl-L-arginyl-L-tryptophylglycyl-L-cysteinyl-L-prolyl-L-prolyl-L-lysyl-L-α-asparagine, trifluoroacetate salt

Synonym: Cyclic [Ac-Cys¹¹, D-Nal¹⁴, Cys¹⁸, Asp-NH₂]-β-MSH(11-22)

Peptide Sequence: Ac-CEHXRWGCPPKD-NH₂ (X = 3-(2-naphthyl)-alanine)

MF: C₇₁H₉₄N₂₀O₁₇S₂ • XCF₃COOH

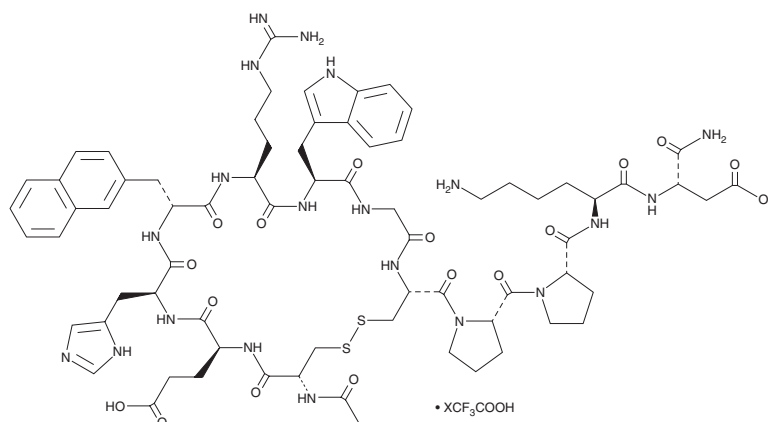
FW: 1,563.8

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

HS-014 (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the HS-014 (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. HS-014 (trifluoroacetate salt) is soluble (≥10 mg/ml) in DMSO.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of HS-014 (trifluoroacetate salt) can be prepared by directly dissolving the solid in aqueous buffers. HS-014 (trifluoroacetate salt) is soluble (≥10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

HS-014 is an antagonist of melanocortin receptor 4 (MC4R; $K_i = 3.16$ nM).¹ It is selective for MC4R over MC1R, MC3R, and MC5R ($K_s = 108, 54.4,$ and 694 nM, respectively). HS-014 inhibits cAMP accumulation induced by α-melanocyte-stimulating hormone (α-MSH) in COS-1 cells expressing MC4R in a concentration-dependent manner. It increases food intake in free-feeding rats when administered intracerebroventricularly at doses of 1 or 3 nmol/animal.² Intracerebroventricular administration of HS-014 (0.008 ng/animal) increases the latency to tail withdrawal in the tail-flick test, as well as inhibits morphine withdrawal-induced decreases in the latency to tail withdrawal in the tail-flick test, in rats.³ It reduces immobility time in the forced swim test, as well as increases time in the open area and reduces the number of droppings in an open field test, in rats when administered at a dose of 100 μg/animal.⁴

References

1. Schiöth, H.B., Mutulis, F., Muceniece, R., et al. *Br. J. Pharmacol.* **124**(1), 75-82 (1998).
2. Kask, A., Rågo, L., Mutulis, F., et al. *Biochem. Biophys. Res. Commun.* **245**(1), 90-93 (1998).
3. Kalange, A.S., Kokare, D.M., Singru, P.S., et al. *Brain Res.* **1181**, 10-20 (2007).
4. Serova, L.I., Laukova, M., Alaluf, L.G., et al. *Behav. Brain Res.* **250**, 139-147 (2013).

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, 08/21/2024

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