

PRODUCT INFORMATION



hMC1R Agonist 1 (acetate)

Item No. 41560

Formal Name: (5R,8S,11S,14R,17S,20R)-17-((1H-imidazol-5-yl)methyl)-8-((1H-indol-3-yl)methyl)-20-((S)-2-acetamidohexanamido)-14-benzyl-11-(3-guanidinopropyl)-7,10,13,16,19-pentaoxo-3,22-dithia-6,9,12,15,18-pentaaza-1(1,3)-benzenacyclotricosaphane-5-carboxamide, acetate

Synonyms: Ac-Nle-[CHfRWC]-NH₂, Ac-Nle-[Cys-His-D-Phe-Arg-Trp-Cys]-NH₂

Peptide Sequence: Ac-XCHfRWC-NH₂
(X=Norleucine)

MF: C₅₄H₇₀N₁₄O₈S₂ • XC₂H₄O₂

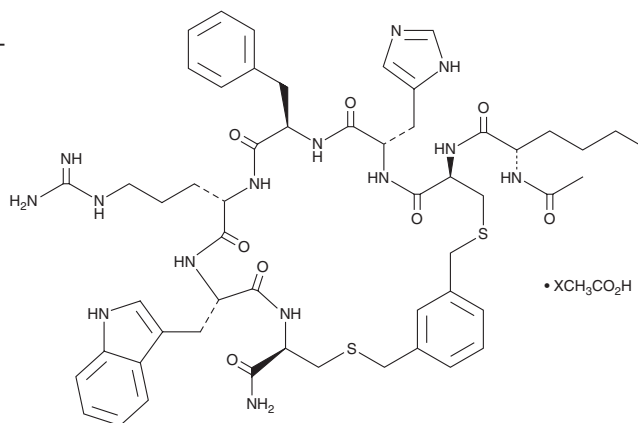
FW: 1,107.4

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

hMC1R agonist 1 (acetate) is supplied as a solid. A stock solution may be made by dissolving the hMC1R agonist 1 (acetate) in the solvent of choice, which should be purged with an inert gas. hMC1R agonist 1 (acetate) is soluble (≥10 mg/ml) in ethanol and DMSO.

Description

hMC1R agonist 1 is a peptide agonist of melanocortin receptor 1 (MC1R) and a derivative of melanotan-II (Item No. 27316).¹ It selectively induces cAMP production in HEK293 cells expressing recombinant human MC1R over those expressing recombinant human MC3R, MC4R, or MC5R (EC₅₀s = 3, 902, 915, and >1,000 nM, respectively).

Reference

1. Tomassi, S., Dimmito, M.P., Cai, M., *et al.* CLIPSing melanotan-II to discover multiple functionally selective hMCR agonists. *J. Med. Chem.* **65**(5), 4007-4017 (2022).

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, 09/11/2024

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