

PRODUCT INFORMATION

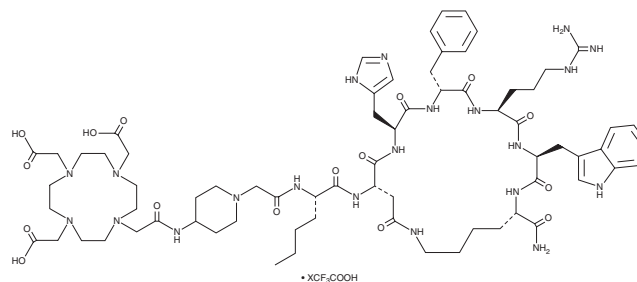


CCZ01048 (trifluoroacetate salt)

Item No. 41459

Formal Name: 2,2',2''-(10-(2-((1-(2-(((S)-1-(((3S,6S,9R,12S,15S,23S)-12-((1H-imidazol-5-yl)methyl)-3-((1H-indol-3-yl)methyl)-9-benzyl-23-carbamoyl-6-(3-guanidinopropyl)-2,5,8,11,14,17-hexaoxo-1,4,7,10,13,18-hexaazacyclotricosan-15-yl)amino)-1-oxohexan-2-yl)amino)-2-oxoethyl)piperidin-4-yl)amino)-2-oxoethyl)-1,4,7,10-tetraazacyclododecane-1,4,7-triyl)triacetic acid, trifluoroacetate salt

Synonym: DOTA-(Pip)-Nle-CycMSH_{hex}
MF: C₇₁H₁₀₅N₂₁O₁₆ • XCF₃COOH
FW: 1,508.7
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

CCZ01048 (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the CCZ01048 (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. CCZ01048 (trifluoroacetate salt) is slightly soluble (0.1-1 mg/ml) in acetonitrile and DMSO.

CCZ01048 (trifluoroacetate salt) is slightly soluble (0.1-1 mg/ml) in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

CCZ01048 is a synthetic peptide composed of a cyclized fragment of α -melanocyte stimulating hormone (α -MSH) conjugated to the bifunctional chelator DOTA.¹ It has been conjugated to radioligands for use in PET imaging of tumors expressing melanocortin receptor 1 (MC1R) and for tissue distribution studies *in vivo*.

Reference

1. Zhang, C., Zhang, Z., Lin, K.S., *et al.* Preclinical melanoma imaging with ⁶⁸Ga-labeled α -melanocyte-stimulating hormone derivatives using pet. *Theranostics* **7**(4), 805-813 (2017).

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, 11/11/2024

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