

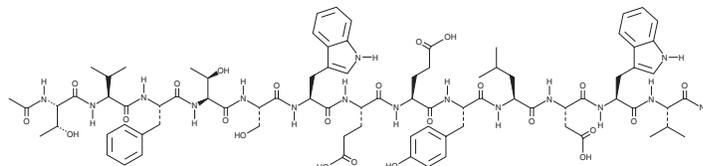
PRODUCT INFORMATION



Pep2-8

Item No. 36923

CAS Registry No.: 1541011-97-5
Formal Name: (3S,6S,9S,12S,15S)-15-((2S,5S,8S,11S,14S,17S)-2-((1H-indol-3-yl)methyl)-11-benzyl-8,17-bis((R)-1-hydroxyethyl)-5-(hydroxymethyl)-14-isopropyl-4,7,10,13,16,19-hexaoxo-3,6,9,12,15,18-hexaazaicosanamido)-3-(((S)-1-(((S)-1-amino-3-methyl-1-oxobutan-2-yl)amino)-3-(1H-indol-3-yl)-1-oxopropan-2-yl)carbamoyl)-12-(2-carboxyethyl)-9-(4-hydroxybenzyl)-6-isobutyl-5,8,11,14-tetraoxo-4,7,10,13-tetrazaooctadecanedioic acid



Synonym: TVFTSWEEYLDWV-NH₂

MF: C₈₃H₁₁₀N₁₆O₂₄

FW: 1,715.9

Purity: ≥95%

UV/Vis.: λ_{max}: 223, 282 nm

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

Pep2-8 is supplied as a solid. A stock solution may be made by dissolving the Pep2-8 in the solvent of choice, which should be purged with an inert gas. Pep2-8 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of Pep2-8 in these solvents is approximately 12 and 5 mg/ml, respectively.

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 Pep2-8 can be prepared by directly dissolving the solid in aqueous buffers. The solubility of Pep2-8 in PBS (pH 7.2) is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Pep2-8 is a peptide inhibitor of proprotein convertase subtilisin kexin type 9 (PCSK9; IC₅₀ = 1.4 μM).¹ It reverses PCSK9-induced decreases in surface LDL receptor levels and restores uptake of LDL in HepG2 cells in a concentration-dependent manner.

Reference

1. Zhang, Y., Eigenbrot, C., Zhou, L., *et al.* Identification of a small peptide that inhibits PCSK9 protein binding to the low density lipoprotein receptor. *J. Biol. Chem.* **289**(2), 942-955 (2014).

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, 12/21/2022

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