

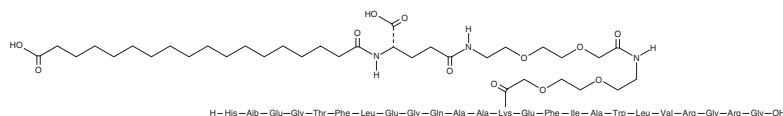
PRODUCT INFORMATION



des-(7-13) Semaglutide

Item No. 44040

Formal Name: (2S,8S,16S,19S,22S,27S,33S,36S,39S,42S,69S)-2-((1H-imidazol-4-yl)methyl)-42-((6S,12S,15S,18S,21S,24S,27S,30S,33S)-21-((1H-indol-3-yl)methyl)-1-amino-30-benzyl-27-((S)-sec-butyl)-12-(3-guanidinopropyl)-33-(hydroxyamino)-6-(2-(hydroxyamino)acetamido)-1-imino-18-isobutyl-15-isopropyl-24-methyl-7,10,13,16,19,22,25,28,31,32-decaoxo-2,8,11,14,17,20,23,26,29-nonaazahexatriacontan-36-amido)-33-(3-amino-3-oxopropyl)-19-benzyl-27-(hydroxyamino)-16-((R)-1-hydroxyethyl)-22-isobutyl-5,5,36,39-tetramethyl-1,6,11,12,15,18,21,24,28,29,32,35,38,41,48,57,66,71-octadeca-50,53,59,62-tetraoxa-3,4,7,14,17,20,23,31,34,37,40,47,56,65,70-pentadecaazaheptaoctacontane-8,69,87-tricarboxylic acid



MF: C₁₅₆H₂₄₆N₃₈O₄₅
FW: 3,373.9
Purity: ≥90%
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

des-(7-13) Semaglutide is supplied as a solid. A stock solution may be made by dissolving the des-(7-13) semaglutide in the solvent of choice, which should be purged with an inert gas. des-(7-13) Semaglutide is slightly soluble (0.1-1 mg/ml) in acetonitrile:water (1:1).

Description

des-(7-13) Semaglutide is a derivative of the glucagon-like peptide 1 receptor (GLP-1R) agonist semaglutide (Item Nos. 40231 | 29969 | 40170) that does not contain amino acids 7-13, which correspond to amino acids 13-19 in GLP-1 and includes two residues important for GLP-1R interaction.^{1,2}

References

1. Adelhors, K., Hedegaard, B.B., Knudsen, L.B., *et al.* Structure-activity studies of glucagon-like peptide-1. *J. Biol. Chem.* **69(9)**, 6275-6278 (1994).
2. Cheang, J.Y. and Moyle, P.M. Glucagon-like peptide-1 (GLP-1)-based therapeutics: Current status and future opportunities beyond type 2 diabetes. *ChemMedChem* **13(7)**, 662-671 (2018).

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, 10/07/2025

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