

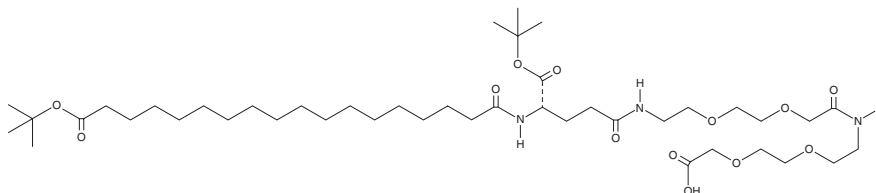
PRODUCT INFORMATION



Semaglutide Side Chain

Item No. 40746

CAS Registry No.: 1118767-16-0
Formal Name: 22S-[(1,1-dimethylethoxy) carbonyl]-10,19,24-trioxo-3,6,12,15-Tetraoxa-9,18,23-triazahentetracontanedioic acid, 41-(1,1-dimethylethyl) ester
MF: $C_{43}H_{79}N_3O_{13}$
FW: 846.1
Purity: $\geq 95\%$
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

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

Description

Semaglutide side chain is a synthetic intermediate.¹ It has been used as an intermediate in the synthesis of the glucagon-like peptide 1 receptor (GLP-1R) agonist semaglutide (Item Nos. 40231 | 29969 | 40170), as well as a lipid side chain to increase the proteolysis-resistance of insulin derivatives for use *in vivo*.^{1,2}

References

1. Guryanov, I., Orlandin, A., De Paola, I., *et al.* Copper(II) lysinate and pseudoproline assistance in the convergent synthesis of the GLP 1 receptor agonists liraglutide and semaglutide. *Org. Process Res. Dev.* **25**(7), 1598-1611 (2021).
2. Kjeldsen, T.B., Hubálek, F., Tagmose, T.M., *et al.* Engineering of orally available, ultralong-acting insulin analogues: Discovery of OI338 and OI320. *J. Med. Chem.* **64**(1), 616-628 (2021).

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, 05/17/2024

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