

# PRODUCT INFORMATION



## Fmoc-L-Lys[Oct-(OtBu)-Glu-(OtBu)-AEEA-AEEA]-OH

Item No. 40707

CAS Registry No.: 1662688-20-1

Formal Name: 3S-carboxy-3OS-[(1,1-dimethylethoxy) carbonyl]-9,18,27,32-tetraoxo-11,14,20,23-tetraoxa-2,8,17,26,31-pentaazonatetracontanedioic acid, 49-(1,1-dimethylethyl) 1-(9H-fluoren-9-ylmethyl) ester

Synonym: Fmoc-Lys(tBuOCO-(CH<sub>2</sub>)<sub>16</sub>-CO-Glu(AEEA-AEEA)-OtBu)-OH

MF: C<sub>64</sub>H<sub>101</sub>N<sub>5</sub>O<sub>16</sub>

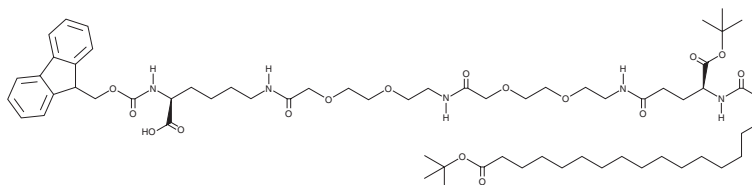
FW: 1,196.5

Purity: ≥95%

Supplied as: A solution in ethanol

Storage: -20°C

Stability: ≥1 year



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

Fmoc-L-Lys[Oct-(OtBu)-Glu-(OtBu)-AEEA-AEEA]-OH is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice, which should be purged with an inert gas. Fmoc-L-Lys[Oct-(OtBu)-Glu-(OtBu)-AEEA-AEEA]-OH is soluble (≥10 mg/ml) in DMSO.

Fmoc-L-Lys[Oct-(OtBu)-Glu-(OtBu)-AEEA-AEEA]-OH is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of Fmoc-L-Lys[Oct-(OtBu)-Glu-(OtBu)-AEEA-AEEA]-OH should be diluted with the aqueous buffer of choice. Fmoc-L-Lys[Oct-(OtBu)-Glu-(OtBu)-AEEA-AEEA]-OH is slightly soluble (0.1-1 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

### Description

Fmoc-L-Lys[Oct-(OtBu)-Glu-(OtBu)-AEEA-AEEA]-OH is a building block in the synthesis of the glucagon-like peptide 1 receptor (GLP-1R) agonist semaglutide (Item Nos. 40231 | 29969 | 40170).<sup>1</sup>

### Reference

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).

#### 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, 05/06/2024

#### 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