

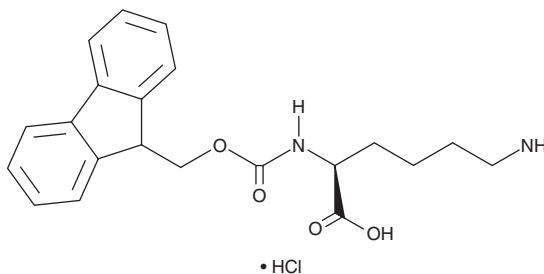
PRODUCT INFORMATION



Fmoc-Lys-OH (hydrochloride)

Item No. 30561

CAS Registry No.: 139262-23-0
Formal Name: N²-[(9H-fluoren-9-ylmethoxy)carbonyl]-L-lysine, monohydrochloride
MF: C₂₁H₂₄N₂O₄ • HCl
FW: 404.9
Purity: ≥98%
UV/Vis.: λ_{max}: 265 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



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

Description

Fmoc-Lys-OH is a building block.^{1,2} It has been used in the synthesis of azido-protected peptides for activity-based protein profiling and click-functional peptide bundles for biopolymer formation.

References

1. Katayama, H., Hojo, H., Ohira, T., *et al.* An efficient peptide ligation using azido-protected peptides via the thioester method. *Tetrahedron Lett.* **49(38)**, 5492-5494 (2008).
2. Wu, D., Sinha, N., Lee, J., *et al.* Polymers with controlled assembly and rigidity made with click-functional peptide bundles. *Nature* **574(7780)**, 658-662 (2019).

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, 05/15/2020

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