

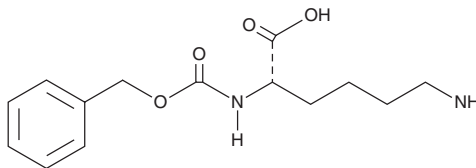
PRODUCT INFORMATION



Z-Lys-OH

Item No. 30460

CAS Registry No.: 2212-75-1
Formal Name: N²-[(phenylmethoxy)carbonyl]-L-lysine
Synonyms: N^α-(Benzyloxycarbonyl)lysine,
α-Carbobenzoxyl-L-lysine
MF: C₁₄H₂₀N₂O₄
FW: 280.3
Purity: ≥98%
Supplied as: A crystalline 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

Z-Lys-OH is supplied as a crystalline solid. Aqueous solutions of Z-Lys-OH can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of Z-Lys-OH in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Z-Lys-OH is an amino acid building block.^{1,2} It has been used in the synthesis of peptidomimetics and polypeptide copolymers.

References

1. Danalev, D., Koleva, M., Ivanova, D.M., *et al.* Synthesis of two peptide mimetics as markers for chemical changes of wool's keratin during skin unhairing process. *Protein Pept. Lett.* **15**(4), 353-355 (2008).
2. Negri, G.E. and Deming, T.J. Protein complexation and pH dependent release using boronic acid containing PEG-polypeptide copolymers. *Macromol. Biosci.* **17**(1), 1600136 (2017).

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, 12/08/2022

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