

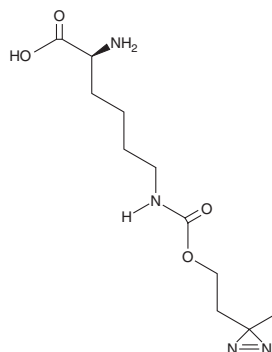
PRODUCT INFORMATION



AbK

Item No. 38895

CAS Registry No.: 1253643-88-7
Formal Name: N⁶-[[2-(3-methyl-3H-diazirin-3-yl)ethoxy]carbonyl]-L-lysine
Synonyms: DiAzKs, H-L-Photo-lysine
MF: C₁₁H₂₀N₄O₄
FW: 272.3
Purity: ≥98%
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

AbK is supplied as a solid. A stock solution may be made by dissolving the AbK in the solvent of choice, which should be purged with an inert gas. AbK is soluble in aqueous acid. It is also slightly soluble in water (warmed). We do not recommend storing the aqueous solution for more than one day.

Description

AbK is an unnatural amino acid¹ It has been incorporated into proteins via a modified aminoacyl-tRNA synthetase and used to cross-link proteins in bacteria or mammalian cells.^{1,2} It has been incorporated into the RNA-binding protein iron regulatory protein 1 (IRP1) and used in the study of RNA-protein interactions *in vitro*.²

References

1. Ai, H.-W., Shen, W., Sagi, A., *et al.* Probing protein-protein interactions with a genetically encoded photo-crosslinking amino acid. *Chembiochem* **12**(12), 1854-1857 (2011).
2. Dziuba, D., Hoffmann, J.-E., Hentze, M.W., *et al.* A genetically encoded diazirine analogue for RNA-protein photo-crosslinking. *Chembiochem* **21**(1-2), 88-93 (2020).

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, 05/28/2024

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