

PRODUCT INFORMATION



DiZPK

Item No. 37191

CAS Registry No.: 1337883-32-5

Formal Name: N⁶-[[[3-(3-methyl-3H-diazirin-3-yl)propyl]amino]carbonyl]-L-lysine

MF: C₁₂H₂₃N₅O₃

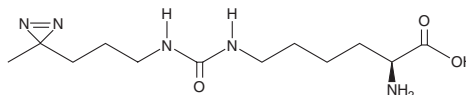
FW: 285.3

Purity: ≥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

DiZPK is supplied as a solid. A stock solution may be made by dissolving the DiZPK in the solvent of choice, which should be purged with an inert gas. DiZPK is slightly soluble in methanol.

Description

DiZPK is a photo cross-linking agent.^{1,2} It contains a diazirine moiety, which, upon UV activation, creates reactive carbenes that form covalent bonds with amino acid side chains and peptide backbones. Incorporation of DiZPK into proteins has been used to map protein-protein interactions in live cells and monitor bacterial defense protein interactions with enteric bacterial pathogens.

References

1. Li, C., Zhang, N., Zhou, J., *et al.* Peptide blocking of PD-1/PD-L1 interaction for cancer immunotherapy. *Cancer Immunol. Res.* **6**(2), 178-188 (2018).
2. Yang, Y., Song, H., and Chen, P.R. Genetically encoded photocrosslinkers for identifying and mapping protein-protein interactions in living cells. *IUBMB Life* **68**(11), 879-886 (2016).

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, 05/16/2023

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