

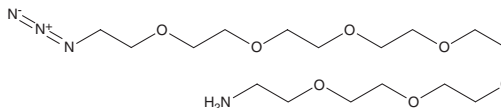
PRODUCT INFORMATION



Azido-dPEG₇-amine

Item No. 30544

CAS Registry No.: 1333154-77-0
Formal Name: 23-azido-3,6,9,12,15,18,21-heptaooxatricosan-1-amine
MF: C₁₆H₃₄N₄O₇
FW: 394.5
Purity: ≥90%
Supplied as: A liquid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

Azido-dPEG₇-amine is supplied as a liquid. A stock solution may be made by dissolving the azido-dPEG₇-amine in the solvent of choice, which should be purged with an inert gas. Azido-dPEG₇-amine is soluble in the organic solvent DMSO at a concentration of approximately 10 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions can be prepared by directly dissolving azido-dPEG₇-amine in aqueous buffers. The solubility of Azido-dPEG₇-amine in PBS, pH 7.2, is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Azido-dPEG₇-amine is a heterobifunctional PEGylated linker.¹ It contains a reactive NH₂ end group, as well as an azide functional group for use in click chemistry reactions. Azido-dPEG₇-amine has been used as a linker in the synthesis of bivalent mammalian target of rapamycin (mTOR) inhibitors.²

References

1. Goswami, L.N., Houston, Z.H., Sarma, S.J., *et al.* Efficient synthesis of diverse heterobifunctionalized clickable oligo(ethylene glycol) linkers: potential applications in bioconjugation and targeted drug delivery. *Org. Biomol. Chem.* **11(7)**, 116-1126 (2012).
2. Rodrik-Outmezguine, V.S., Okaniwa, M., Yao, Z., *et al.* Overcoming mTOR resistance mutations with a new-generation mTOR inhibitor. *Nature* **534(7606)**, 272-276 (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, 05/11/2020

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