

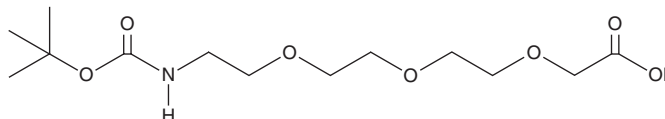
PRODUCT INFORMATION



Boc-NH-PEG₃-CH₂COOH

Item No. 30536

CAS Registry No.: 462100-06-7
Formal Name: 5,8,11-trioxa-2-azatridecanedioic acid, 1-(1,1-dimethylethyl) ester
MF: C₁₃H₂₅NO₇
FW: 307.3
Purity: ≥95%
Supplied as: A neat oil
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

Boc-NH-PEG₃-CH₂COOH is supplied as a neat oil. A stock solution may be made by dissolving the Boc-NH-PEG₃-CH₂COOH in the solvent of choice, which should be purged with an inert gas. Boc-NH-PEG₃-CH₂COOH is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of Boc-NH-PEG₃-CH₂COOH in these solvents is approximately 30, 1, and 5 mg/ml, respectively.

Description

Boc-NH-PEG₃-CH₂COOH is a building block.^{1,2} It has been used as a linker in the synthesis of proteolysis-targeting chimeras (PROTACs) targeting MEK.¹ It has also been used as a spacer in the synthesis of radiolabeled probes for dual labeling of tumor integrin αVβ3 and the gastrin-releasing peptide receptor (GRPR) in tumor tissue *in vivo*.²

References

1. Vollmer, S., Cunoosamy, D., Lv, H., *et al.* Design, synthesis, and biological evaluation of MEK PROTACs. *J. Med. Chem.* **63**(1), 157-162 (2020).
2. Liu, Z., Yan, Y., Chin, F.T., *et al.* Dual integrin and gastrin-releasing peptide receptor targeted tumor imaging using ¹⁸F-labeled PEGylated RGD-bombesin heterodimer ¹⁸F-FB-PEG₃-Glu-RGD-BBN. *J. Med. Chem.* **52**(2), 425-432 (2009).

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, 05/19/2020

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