

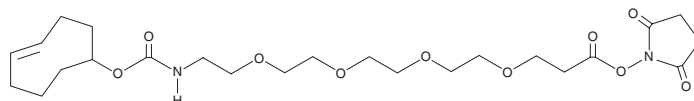
PRODUCT INFORMATION



TCO-PEG₄-NHS ester

Item No. 44735

CAS Registry No.: 1621096-79-4
Formal Name: 5,8,11,14-tetraoxa-2-azaheptadecanedioic acid, 1-(4E)-4-cycloocten-1-yl 17-(2,5-dioxo-1-pyrrolidinyl) ester
Synonym: *trans*-Cyclooctene-PEG₄-N-hydroxysuccinimide
MF: C₂₄H₃₈N₂O₁₀
FW: 514.6
Purity: ≥95%
Supplied as: A 50 mg/ml solution in acetonitrile
Storage: -20°C
Stability: ≥4 years



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

Description

TCO-PEG₄-NHS ester is a click chemistry building block.¹ It has been used as a starting material in the synthesis of HELIOS-based fluorescent probes and radiotracers for *in vivo* imaging of bombesin receptor subtype 2-expressing prostate cancer mouse xenograft models.^{1,2}

References

1. Meimetis, L.G., Carlson, J.C.T., Giedt, R.J., *et al.* Ultrafluorogenic coumarin-tetrazine probes for real-time biological imaging. *Angew Chem. Int. Ed. Engl.* **53**(29), 7531-7534 (2014).
2. Handula, M., Verhoeven, M., Chen, K.T., *et al.* Towards complete tumor resection: Novel dual-modality probes for improved image-guided surgery of GRPR-expressing prostate cancer. *Pharmaceutics* **14**(1), 195 (2022).

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, 05/12/2026

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