

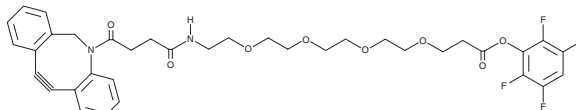
PRODUCT INFORMATION



DBCO-PEG₄-TFP ester

Item No. 45967

Synonym: Dibenzocyclooctyne-PEG₄-tetrafluorophenyl ester
MF: C₃₆H₃₆F₄N₂O₈
FW: 700.7
Purity: ≥95%
Supplied as: A 10 mg/ml solution in ethanol
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

DBCO-PEG₄-TFP ester is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. DBCO-PEG₄-TFP ester is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in methanol.

Description

DBCO-PEG₄-TFP ester is a labeling reagent containing an amine-reactive tetrafluorophenyl (TFP) ester moiety and an azido-reactive dibenzocyclooctyne (DBCO) copper-free click chemistry moiety connected via a PEG linker. It has been used to conjugate antibodies to metal-chelating polymers for mass cytometry.¹

Reference

1. Zhang, Y., Liu, P., Majonis, D., *et al.* Polymeric dipicolylamine based mass tags for mass cytometry. *Chem. Sci.* **13**(11), 3233-3243 (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, 07/20/2026

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