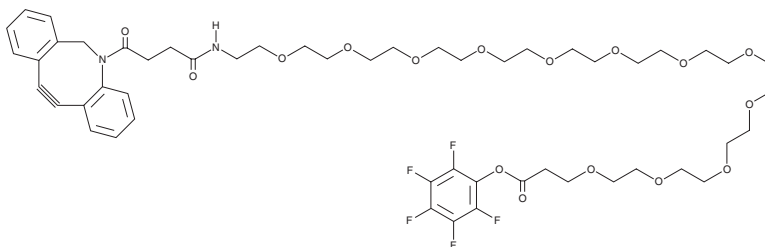


PRODUCT INFORMATION



DBCO-PEG₁₂-PFP ester Item No. 45971

Synonym: Dibenzocyclooctyne-PEG₁₂-pentafluorophenyl ester
MF: C₅₂H₆₇F₅N₂O₁₆
FW: 1,071.1
Purity: ≥90%
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-PEG₁₂-PFP ester is supplied as a solution in ethanol. A stock solution may be made by dissolving the DBCO-PEG₁₂-PFP ester in the solvent of choice, which should be purged with an inert gas. DBCO-PEG₁₂-PFP ester is slightly soluble (0.1-1 mg/ml) in methanol.

Description

DBCO-PEG₁₂-PFP ester is a labeling reagent containing an amine-reactive pentafluorophenyl (PFP) ester moiety and an azido-reactive dibenzocyclooctyne (DBCO) copper-free click chemistry moiety connected via a PEG linker.

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, 07/14/2026

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