

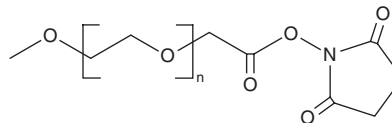
PRODUCT INFORMATION



mPEG(30000)-NHS ester

Item No. 39743

Formal Name: 2,5-dioxopyrrolidin-1-yl
2-(2-methoxyethoxy)acetate
Synonyms: Methoxy Polyethylene Glycol-30000-
NHS ester, MPEG₃₀₀₀₀-NHS ester
MF: (C₂H₄O)_nC₇H₉NO₅
Purity: ≥90%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

mPEG(30000)-NHS ester is supplied as a solid. A stock solution may be made by dissolving the mPEG(30000)-NHS ester in water. We do not recommend storing the aqueous solution for more than one day.

Description

mPEG(30000)-NHS ester is a PEGylated and esterified form of N-hydroxysuccinimide (NHS).¹

Reference

1. Meziadi, A., Zuberi, N., de Haan, H.W., *et al.* Overcoming PEG—protein mutual repulsion to improve the efficiency of PEGylation. *Biomacromolecules* **23(11)**, 4948-4956 (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, 04/11/2025

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