

PRODUCT INFORMATION



Biotin-mPEG(2000)

Item No. 46013

CAS Registry No.: 767356-08-1

Formal Name: α -[2-[[5-[(3aS,4S,6aR)-hexahydro-2-oxo-1H-thieno[3,4-d]imidazol-4-yl]-1-oxopentyl]amino]ethyl]- ω -methoxy-poly(oxy-1,2-ethanediyl)

Synonym: PEG_{2k}-biotin

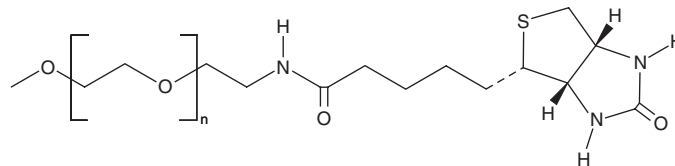
MF: (C₂H₄O)_nC₁₃H₂₃N₃O₃S

Purity: $\geq 90\%$

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

Biotin-mPEG(2000) is supplied as a solid. A stock solution may be made by dissolving the biotin-mPEG(2000) in the solvent of choice, which should be purged with an inert gas. Biotin-mPEG(2000) is soluble (≥ 10 mg/ml) in DMSO and sparingly soluble (1-10 mg/ml) in ethanol.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of biotin-mPEG(2000) can be prepared by directly dissolving the solid in aqueous buffers. Biotin-mPEG(2000) is soluble (≥ 10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

Biotin-mPEG(2000) is a surface modification agent.¹ It has been used to modify the surface of mesenchymal stem cells (MSCs) to prevent their adhesion to primary mouse aortic endothelial cells.

Reference

1. Takayama, Y., Kusamori, K., Katsurada, Y., *et al.* Efficient delivery of mesenchymal stem/stromal cells to injured liver by surface PEGylation. *Stem Cell Res. Ther.* **14**(1), 216 (2023).

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, 06/12/2026

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