

PRODUCT INFORMATION



Biotin-PEG₂-Maleimide

Item No. 44569

CAS Registry No.: 305372-39-8
Formal Name: (3a*S*,4*S*,6a*R*)-*N*-[2-[2-[2-[[3-(2,5-dihydro-2,5-dioxo-1*H*-pyrrol-1-yl)-1-oxopropyl]amino]ethoxy]ethoxy]ethyl]hexahydro-2-oxo-1*H*-thieno[3,4-*d*]imidazole-4-pentanamide

Synonym: Biotin-PEG₂-Mal

MF: C₂₃H₃₅N₅O₇S

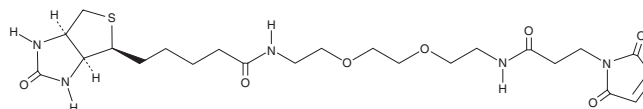
FW: 525.6

Purity: ≥95%

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-PEG₂-maleimide is supplied as a solid. A stock solution may be made by dissolving the biotin-PEG₂-maleimide in the solvent of choice, which should be purged with an inert gas. Biotin-PEG₂-maleimide is soluble (≥10 mg/ml) in DMSO.

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-PEG₂-maleimide can be prepared by directly dissolving the solid in aqueous buffers.

Biotin-PEG₂-maleimide 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-PEG₂-maleimide is a thiol- and alkene-reactive biotinylation reagent.^{1,2} It has been used in the biotinylation of proteins and oligonucleotides.

References

1. Phan, J.C., Nehilla, B.J., Srinivasan, S., *et al.* Human immunodeficiency virus (HIV) separation and enrichment via the combination of antiviral lectin recognition and a thermoresponsive reagent system. *Pharm. Res.* **33(10)**, 2411-2420 (2016).
2. Okamoto, A., Taiji, T., Tainaka, K., *et al.* Oligonucleotides containing 7-vinyl-7-deazaguanine as a facile strategy for expanding the functional diversity of DNA. *Bioorg. Med. Chem. Lett.* **12(15)**, 1895-1896 (2002).

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, 12/17/2025

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