

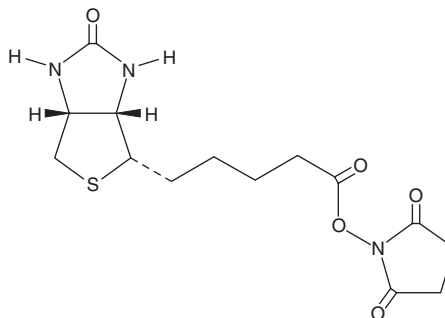
# Product Information



## Biotin-NHS

Catalog No. 13315

**CAS Registry No.:** 35013-72-0  
**Formal Name:** 6aR-hexahydro-2-oxo-2,5-dioxo-1-pyrrolidinyl ester-1H-thieno[3aS,4-d]imidazole-4S-pentanoic acid  
**Synonyms:** Biotin N-Hydroxysuccinimide Ester, NSC 345668  
**MF:** C<sub>14</sub>H<sub>19</sub>N<sub>3</sub>O<sub>5</sub>S  
**FW:** 341.4  
**Purity:** ≥95%  
**Stability:** ≥2 years at -20°C  
**Supplied as:** A crystalline solid



For long term storage, we suggest that biotin-NHS be stored as supplied at -20°C. It should be stable for at least two years.

Biotin-NHS is supplied as a crystalline solid. A stock solution may be made by dissolving the biotin-NHS in the solvent of choice. Biotin-NHS is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of biotin-NHS in these solvents is approximately 20 mg/ml.

Biotin-NHS is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, biotin-NHS should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Biotin-NHS has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

N-Hydroxysuccinimide (NHS) activates carboxylic acid groups (on biotin) to facilitate coupling reactions. Biotin-NHS is a compound used to attach biotin to primary amines under alkaline conditions (pH-8-9). For example, lysines on the surface of proteins, including antibodies, are ideal targets for biotinylation with this compound.

### Laboratory Procedures

Avoid performing biotinylation reactions in buffers containing primary amines (such as Tris). The primary amines in the buffers will compete with the protein for biotinylation. Prior to biotinylation, exchange the buffer to one which does not contain primary amines (such as potassium phosphate). This can be accomplished by dialysis or by running the protein solution over a desalting column. The optimal level of biotinylation will be protein dependent. The procedure below uses a twenty-fold molar excess of NHS-biotin relative to protein, which is appropriate for many proteins. If you are unsure if this level of biotinylation is appropriate for your protein, it may be advisable to set up a number of small biotinylation reactions, increasing and decreasing the ratio of biotinylation reagent to protein by four-fold (i.e., using molar excesses of NHS-biotin to protein of five-fold, twenty-fold and eighty-fold).

1. Dilute protein in 100 mM potassium phosphate, pH 7.4 to a final concentration of 0.5-10 mg/ml.
2. Just prior to use, prepare a 20 mM stock solution of NHS-biotin in dimethyl formamide or DMSO. This solution may be stored at -20°C for up to two weeks; long term storage is not recommended.
3. Add a twenty-fold molar excess of NHS-biotin to the protein.
4. Incubate one hour at room temperature, or overnight at 4°C.
5. Unreacted NHS-biotin may be quenched by the addition of an excess of 1 M Tris, pH 8.0. Alternatively, unreacted NHS-biotin may be removed by dialysis, or purifying with a desalting column.
6. The biotinylated protein is now ready for use.

### Related Products

Biotin-XX hydrazide - Cat. No. 13304 • Biotin-X-NHS - Cat. No. 13316

**WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.**

### MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent via email to your institution.

### WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery.**

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees. Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy** located on our website and in our catalog.

Copyright Cayman Chemical Company, 07/09/2010

## Cayman Chemical

**Mailing address**  
1180 E. Ellsworth Road  
Ann Arbor, MI  
48108 USA

**Phone**  
(800) 364-9897  
(734) 971-3335

**Fax**  
(734) 971-3640

**E-Mail**  
custserv@caymanchem.com

**Web**  
www.caymanchem.com