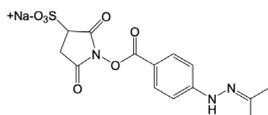


Sulfo S-HyNic Linker (Water Soluble)

C₁₄H₁₄N₃NaO₇S; Mol. Wt.: 391.33

Cat. No.: S-1011-010

Storage: Desiccated: -15° to -25°C



Introduction

SoluLINK® bioconjugation technology is based on the formation of a stable covalent bond that has a UV-traceable signal to indicate the real-time formation of conjugate. This bond is a bis-aryl hydrazone formed from an aromatic hydrazine and an aromatic aldehyde. **Sulfo S-HyNic** (sulfosuccinimidyl 6-hydrazinonicotinate acetone hydrazone) is a water-soluble version of S-HyNic which is used to incorporate aromatic hydrazine linkers on biomolecules and surfaces. Sulfo S-HyNic is an amine-reactive linker that directly converts amino groups to HyNic groups. **S-4FB** (succinimidyl 4-formylbenzoate) is used to convert amino groups to aromatic aldehydes (4-formylbenzamide, or 4FB groups). Addition of a HyNic-modified biomolecule to a 4FB-modified biomolecule or surface leads directly to the formation of the conjugate (Figure 1). The conjugate bond is stable to 92°C and pH 2.0-10.0. The recommended pH for biomolecule modification is 8.0, and conjugation is 6.0. Unlike thiol-based conjugation protocols where reducing reagents are required that can compromise the activity of proteins by cleaving disulfide bonds, the HyNic-4FB conjugation couple leaves disulfide bonds intact. No oxidants, reductants or metals are required in the preparation of conjugate.

Further enhancing the many advantages of the HyNic/4FB conjugation couple is the discovery *Dirksen et al.* that showed aniline catalyzes the formation of this Schiff's base. This is especially effective for large biomolecule conjugations. In the case of antibody-protein conjugations the addition of 10 mM **TurboLINK Catalyst Buffer (10X)** (aniline) to the reaction converts >95% of the antibody to conjugate in ~2 hours using 1 – 2 mole equivalents of the second protein.

The HyNic-4FB conjugation couple is chromophoric - the conjugate bond absorbs at 354 nm and has a molar extinction coefficient of 29,000 L/(mol*cm). This allows (1) real time spectrophotometric monitoring of a conjugate reaction, (2) ability to 'visualize' the conjugate during chromatographic purification using a UV or photodiode array detector and (3) quantification of conjugation. Furthermore, the level of incorporation of HyNic groups can be quantified colorimetrically as reaction with **2-Sulfobenzaldehyde** yields a chromophoric product that absorbs at 350 nm with a molar extinction coefficient of 28,500 L/(mol*cm).

Sulfo S-HyNic is a water-soluble sulfo NHS ester linker that converts amines on biomolecules and surfaces to HyNic groups. Sulfo-S-HyNic is recommended for modification of any amino surfaces such as beads and quantum dots.

Links to procedures and calculators are provided in this user guide. (Also see [Application Notes](#)).

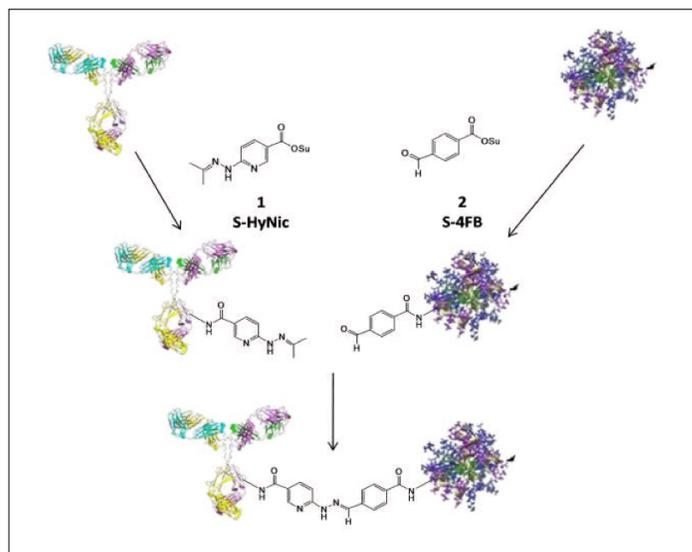


Figure 1: Schematic representation of SoluLINK bioconjugation chemistry where an antibody is modified with S-HyNic to incorporate HyNic groups and a second protein is modified with S-4FB to incorporate 4FB groups. Conjugate is formed directly by simply mixing the HyNic-modified antibody with the 4FB-modified protein.

Additional materials required

Reagents

Zeba™ Desalting Columns
Modification Buffer (10X)
Conjugation Buffer (10X)
Anhydrous DMF
2-Sulfobenzaldehyde

Equipment

Variable-speed bench-top microcentrifuge
Spectrophotometer or Plate Reader
1.5 mL microcentrifuge tubes

Modification Procedure

A. Desalting

- Desalt/buffer exchange the protein into 1X Modification Buffer (100 mM sodium phosphate, 150 mM sodium chloride, pH 8.0). If needed, refer to the [Protein Desalting Protocol](#) or [Oligonucleotide Desalting Protocol](#).

Notes:

- It is necessary to remove all free amine-containing contaminants, e.g., tris or glycine, from the protein before modification.
- High-level buffering capacity, i.e. 100 mM phosphate, is necessary for successful modification.
- For desalting proteins, Zeba Desalting Columns are recommended; for oligonucleotides, Sartorius Vivaspin diafiltration units. Refer to desalting protocol for either apparatus.

Continued on next page.

S-1011-010, LBL-02186.Rev.00

Continued from page 1.

B. Determine the concentration of the desalted protein

1. Determine the concentration of the protein to be modified using a [Bradford assay](#) or [BCA assay](#) or oligonucleotide by using the [Oligonucleotide Concentration Determination Protocol](#). Alternatively the A280 can be used if the protein extinction coefficient is known (E1%).
2. Adjust the concentration to 1.0 – 2.5 mg/mL in 1X Modification Buffer, pH 8.0, if necessary.

C. Prepare Sulfo S-HyNic/DMF stock solution

1. Prepare a stock solution of Sulfo-S-HyNic in anhydrous DMF or aqueous buffer by dissolving 2 – 4 mg of Sulfo-S-HyNic in 100 μ L anhydrous DMF or in aqueous buffer.

Note:

a) The Sulfo-S-HyNic/DMF stock solution is stable for 2 weeks if prepared with anhydrous DMF and stored desiccated.

b) Stock solution prepared in buffer must be used immediately.

D. Modification of a protein

1. Using Table 1 as a guide, add the required volume of Sulfo S-HyNic solution to the protein solution. Refer to the [Protein Modification Calculator](#), or [Amino-Oligonucleotide Modification Calculator](#), if needed.

IgG Concentration (mg/mL)	Sulfo S-HyNic Mole Equivalents Added	Determined Ratio of HyNic/Protein
1.0	20	5.5
	30	8.2
4.0	15	4.7
	20	6.4
	25	7.8

Table 1: The number of HyNic groups incorporated on an antibody is dependent on the number of mole equivalents sulfo-S-HyNic added and the protein concentration. This table can be used as a general guide for modification of any protein with a succinimidyl-based modification reagent.

2. Allow the reaction to incubate at room temperature for 2.0 hours.

E. Desalting procedure

1. Desalt/buffer exchange the protein into 1X Conjugation Buffer (100 mM sodium phosphate, 150 mM sodium chloride, pH 6.0). If needed, refer to the [Protein Desalting Protocol](#) or [Oligonucleotide Desalting Protocol](#).

F. Quantifying modification level

1. The molar substitution ratio (MSR) can be determined using a colorimetric reaction outlined in Figure 2. Addition of 2-sulfobenzaldehyde to a HyNic-modified biomolecule yields a bis-aryl hydrazone that absorbs at 350 nm. Refer to the [HyNic-Protein MSR Calculator](#) as well as the protocol that is appropriate for your lab equipment: [HyNic Protein MSR Instructions](#).

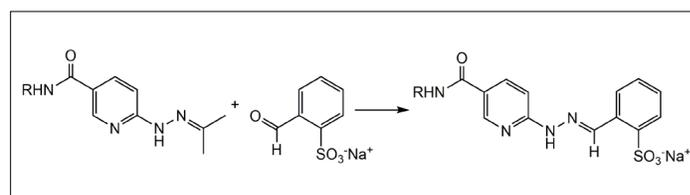


Figure 2: Colorimetric reaction used to quantify number of HyNic moieties on a biomolecule.

2. The biomolecule is now HyNic-modified and ready for conjugation to 4FB-modified biomolecules and surfaces.

Application Notes

[Performing a Bradford assay](#)

[Performing a BCA protein assay](#)

[Troubleshooting Guide](#)