

## Introduction to Solulink Bioconjugation Technology

Solulink's core technology is based on the formation of a stable aromatic bond that has a UV-traceable signal to indicate the real-time formation of the conjugate. S-HyNic **1** (succinimidyl 6-hydrazinonicotinate acetone hydrazone, SANH) is used to incorporate aromatic hydrazine linkers on biomolecules. S-HyNic is an amino-reactive reagent that directly converts amino groups on biomolecules and surfaces to HyNic groups. S-4FB **2** (succinimidyl 4-formylbenzoate, SFB) is used to convert amino groups to aromatic aldehydes (4-formylbenzamide (4FB) groups). Addition of a HyNic-modified biomolecule to a 4FB-modified biomolecule or surface directly leads 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 antibody 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 by Dirksen *et al.*<sup>1</sup> that showed that 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 (aniline) to the reaction mixture converts >95% of the antibody to conjugate in ~2 hours using 1-2 mole equivalents of second protein.

The HyNic-4FB conjugation couple is chromophoric- the conjugate bond absorbs at 354 nm and has a molar extinction coefficient of 29000. 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 4FB groups can be quantified colorimetrically as reaction with 2-Hydrazinopyridine (Solulink catalog# # S-2002) yields a chromophoric product that absorbs at A354 with a molar extinction coefficient of 28,500 (Figure 2). Links to procedures and calculators to guide users through this process are given in the procedures below.

S-SS-4FB is an analog of S-4FB with a disulfide cleavable moiety in the linker if required for intracellular cleavable of the conjugate.

## Methods

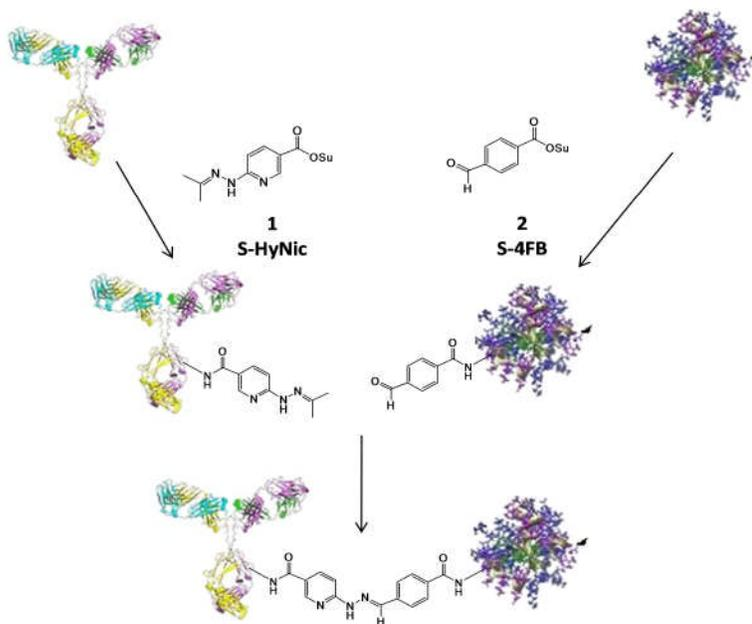
### Additional Materials Required

#### Reagents

Diafiltration spin columns	S-4004
Modification Buffer	S-4003
Conjugation Buffer	S-4002
Anhydrous DMF	S-4001

#### Equipment

Variable-speed bench-top microcentrifuge
Spectrophotometer or Plate Reader
1.5 mL microcentrifuge tubes
2-Sulfobenzaldehyde (S-2005)



**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 proteins.

**Note:** This protocol and all documents linked below can be downloaded from the appropriate category in the Solulink Library at <http://www.solulink.com/library>.

## Modification Procedure

### A. Desalting procedure

- Desalt/buffer exchange the protein or oligonucleotide into 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](#), if needed.

#### Notes:

- It is necessary to remove all free amine-containing contaminants, e.g. tris, glycine, from the protein solution before modification.
- High-level buffering capacity, i.e. 100 mM phosphate, is necessary for successful modification.
- For desalting proteins Solulink recommends Pierce Zeba Desalt Spin columns (# 89882); for oligonucleotides, Sartorius VivaSpin diafiltration units (#VS0112). Refer to desalting protocol for either apparatus.

### B. Determine the concentration of the protein

- Determine the concentration of the protein to be modified using a [Bradford assay](#) (BioRad, #500-0006) or the [BCA assay](#) (ThermoScientific, #23223). Alternatively the A280 can be used if the protein extinction coefficient is known (E1%).
- Adjust the concentration to 1-2.5 mg/mL in Modification Buffer pH 8.0, if necessary

### C. Prepare a S-SS-4FB/DMF stock solution

- Prepare a stock solution of S-SS-4FB in anhydrous DMF (or DMSO).

**Note:** The S-SS-4FB/DMF stock solution is stable for 2 weeks if prepared with anhydrous DMF (Solulink Cat. No. S-4001) and stored desiccated. We recommend dissolving 1 mg of S-SS-4FB in 100  $\mu$ L DMF and 10 mg of S-SS-4FB in 500  $\mu$ L DMF.

### D. Modification of protein or oligonucleotides

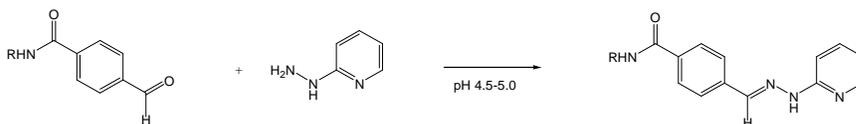
- Using Table 1 as a guide, add the requisite volume of S-SS-4FB/DMF to the protein solution; refer to the [Protein Modification With An NHS Ester Calculator](#) or [Amino-Oligonucleotide Modification With An NHS Ester Calculator](#), if needed.
  - Note:** be sure to use the correct values for S-SS-4FB in the **Reagent Information** section of the calculator
- Allow reaction to incubate at room temperature for 2.0 hours.

### E. Desalting procedure

- Desalt/buffer exchange the protein or oligonucleotide into 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](#), if needed.

### F. Quantifying modification level

- The molar substitution ratio (MSR) can be determined using a colorimetric reaction outlined in Figure 2. Addition of 2-hydrazinopyridine to a 4FB-modified biomolecules yields a compound that absorbs at 350 nm. Refer to the [4FB-Protein Colorimetric MSR Calculator](#) or the [4FB-Oligonucleotide Colorimetric MSR Calculator](#) as well as the protocol that is appropriate for your lab equipment: [4FB Colorimetric MSR Assay Protein Nanodrop Method Protocol](#), [4FB Colorimetric MSR Assay Protein Spectrophotometer Method Protocol](#), [4FB Colorimetric MSR Assay Oligonucleotide Nanodrop Method Protocol](#), or [4FB Colorimetric MSR Assay Oligonucleotide Spectrophotometer Method Protocol](#).



**Figure 2:** Colorimetric reaction used to quantify number of 4FB linkers on a biomolecule

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

## Troubleshooting

Problem	Possible Cause	Solution
Protein was not modified or poorly modified.	Protein has been contaminated with amine containing compounds  The concentration of the protein was too low	Desalt the protein more thoroughly with a new Zeba Spin column or VivaSpin diafiltration apparatus  Increase the concentration of the protein to >2.0 mg/mL
S-SS-4-FB was hydrolyzed	Wet or poor quality DMF/DMSO hydrolyzed the NHS ester	Use a good quality anhydrous DMF/DMSO to solubilize S-SS-4-FB.

## Stability

The SS-4FB linkers incorporated on biomolecules are stable at 4°C for >30 days.

IgG concentration (mg/mL)	S-SS-4FB molar equivalents added	Determined ratio of 4FB/protein
1.0	5	2.38
	10	4.73
	15	6.20
2.5	5	3.08
	10	6.58
5.0	5	3.74
	10	6.80

**Table 1:** The number of 4FB groups incorporated on an antibody is dependent on the number of mole equivalents S-SS-4FB 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.

## Related SoluLink Products

S-4004	VivaSpin diafiltration device	S-2002	2-Hydrazinopyridine	S-2006	TurboLink Catalyst Buffer
S-1004	S-4FB Linker	S-4001	Anhydrous DMF		
S-4003	Modification Buffer	S-4002	Conjugation Buffer		

## References

1. Dirksen, A., et al., Nucleophilic catalysis of hydrazone formation and transimination: implications for dynamic covalent chemistry. *J Am Chem Soc*, 2006. 128(49): p. 15602-3.
2. Ryan C. Bailey, Gabriel A. Kwong, Caius G. Radu, Owen N. Witte, and James R. Heath, DNA-Encoded Antibody Libraries: A Unified Platform for Multiplexed Cell Sorting and Detection of Genes and Proteins, *J. Amer. Chem. Soc.* 2007, **129**, 1959-1967.

The products offered here are for research use only. Any commercial application will require a license from Solulink. The Solulink Conjugation System is patented and has multiple patents pending. Please contact Solulink for information regarding licensing information. Solulink products and methods may be covered by one or more of the following United States patents Nos. 6,686,461, 6,800,728, 7,102,024, 7,173,125, 7,462,689 and other pending patent applications. Information in this manual is subject to change without notice and does not constitute a commitment on the part of Solulink, Inc. It is supplied on an "as is" basis without any warranty of any kind, either explicit or implied. Information may be changed or updated in this manual at any time. This document may not be copied, transferred, reproduced, disclosed, or duplicated, in whole or in part, without the prior written consent of Solulink, Inc. This documentation is proprietary information and protected by the copyright laws of the United States and international treaties. The manufacturer of this documentation is Solulink, Inc.