

Modify and label oligonucleotide 5' phosphate groups

TR0030.5

Introduction

The 5' phosphate group of oligonucleotides, DNA and RNA can be conjugated to primary amine-containing molecules using the carbodiimide crosslinker EDC (Part No. 22980) and imidazole. The example protocol presented in this Tech Tip describes the amine-modification of an oligonucleotide with an excess of ethylenediamine, as illustrated in Figure 1. Depending on the amine-containing molecules used, the crosslinking strategy can be adapted in a number of ways to directly or indirectly modify, label or conjugate an oligonucleotide (Table 1).



Figure 1. Amine-modification of oligonucleotide using ethylenediamine, EDC and imidazole.

Conjugation Goal	Reagent (Thermo Scientific Part No.) and Strategy
Biotin-label the oligonucleotide	Use Biotin Hydrazide (21339) or Biotin-LC-Hydrazide (21340) instead of ethylenediamine in the default reaction.
Create a sulfhydryl-reactive oligonucleotide	Use EMCH (22106), KMUH (22111) or MPBH (22305) instead of ethylenediame in the default reaction. Use PDPH (22301) instead of ethylenediamine to obtain a sulfhydryl crosslink that is reversible. This strategy is useful for preparing conjugates with sulfhydryl-containing proteins and other molecules.
Create a sulfhydryl group on oligonucleotide	Use cystamine (H ₂ N-CH ₂ -CH ₂ -S-S-CH ₂ -CH ₂ -NH ₂) instead of ethylenediamine in the default reaction, and then reduce the disulfide bond with DTT or similar reagent. This strategy is useful for preparing oligo-enzyme conjugates (by reaction to maleimide-activated enzymes, such as Part No. 31485) for assays and blotting procedures.
Immobilize oligonucleotide to a beaded affinity support	Use UltraLink Hydrazide (53149) instead of ethylenediamine in the default reaction. This strategy is useful for preparing supports for affinity purification of binding partners.
Fluorophore-label the oligonucleotide	Amine-modify the oligo with ethylenediamine using the default procedure, then react the amino group with the desired amine-reactive fluorophore (e.g., Fluorescein, 46410 or 46425; DyLight [™] 550, 62262; or DyLight 650, 62265)

Table 1. Possible applications for EDC/imidazole-mediated modification of oligonucleotides.



Materials Required

- EDC (1-ethyl-3-[3-dimethylaminopropyl]carbodiimide hydrochloride), Part No. 22980
- Ethylenediamine (Part No. 23031) or other amine- or hydrazide-containing molecule
- 0.1 M Imidazole, pH 6
- Reaction Buffer, such as phosphate buffered saline (PBS) with EDTA: 10 mM sodium phosphate, 0.15M NaCl, 10mM EDTA, pH 7.2. Avoid using PBS with > 10mM phosphate, which will interfere with the intended reaction. Other amine-free and carboxylate-free buffers may be substituted, but avoid Tris, which contains a primary amine that will quench the reaction.
- 7.5-15nmol (~60-120µg) oligonucleotide or double-stranded DNA or RNA dissolved in ~10µL Reaction Buffer.

Procedure for Modifying 5' Phosphate Groups

- 1. Dissolve ethylenediamine (or alternative) to a final concentration of 0.25M in 10µL of 0.1M imidazole.
- 2. Weigh 1.25mg (6.52µmol) of EDC into a microcentrifuge tube.
- 3. Add 7.5 μ L of the prepared oligonucleotide to the tube containing the EDC and immediately add 5 μ L of the ethylenediamine/imidazole solution.
- 4. Vortex tube until contents are completely dissolved, and then briefly centrifuge the tube to gather contents.
- 5. Add an additional 20µL of 0.1M imidazole, pH 6.
- 6. Incubate reaction at one of the following conditions:

Temperature	Time
50°C	30 minutes to 2 hours
37°C	1 hour to overnight
Room Temperature	2 hours to overnight

 Remove non-reacted EDC and its by-products and imidazole by dialysis (e.g., Slide-A-Lyzer[®] MINI Dialysis Units) or spin desalting column (Zeba[™] Spin Desalting Column) using 10mM sodium phosphate, 0.15M NaCl, 10mM EDTA, pH 7.2, or other suitable buffer.

Note: The amine-modified oligonucleotide may be stored frozen for up to one year or used immediately for other conjugation reactions. If heterobifunctional hydrazide compounds were used, purify and store the conjugate in a manner suitable for stability of the second reactive group.

Additional Information

The following reference (Part No. 20036) provides further discussion of EDC-mediated oligonucleotide modification and other oligonucleotide crosslinking techniques, as well as citations of original literature:

Hermanson, G.T. (2008). Bioconjugate Techniques. 2nd Edition. Academic Press, San Diego.

© 2011 Thermo Fisher Scientific Inc. All rights reserved. Unless otherwise indicated, all trademarks are property of Thermo Fisher Scientific Inc. and its subsidiaries. Printed in the USA.