Labeling Azide-Modified Biomolecules with Fluorescent Dye Alkynes

Labeling Oligonucleotides with Dye Alkynes

- 1. Prepare the following stock solutions:
 - 200 mM THPTA [tris(3-hydroxypropyltriazolylmethyl)amine)] in water
 - 100 mM CuSO4 in water
 - Azide-modified oligo in water (as concentrated as possible, e.g., >10 mg/ml)
 - 100 mM sodium ascorbate in water
 - 10 mM dye alkyne in DMSO or water (see our website for recommended solvent)
- 2. Mix and vortex well CuSO4 with THPTA in a 1:2 ratio for several minutes before the reaction. This working solution is stable for several weeks when frozen.
- **3**. To the azide-modified oligo solution, add an excess of dye alkyne (2-5 equivalents by molar ratio).
- 4. Add 5 equivalents of THPTA/CuSO4 working solution (from Step 1)
- 5. Add 10-30 equivalents of sodium ascorbate.
- 6. Stir, vortex or shake the reaction mixture at room temperature for 30-60 minutes.
- 7. Ethanol-precipitate the oligo or purify it by your desired method (e.g., HPLC).

Labeling Peptides with Dye Alkynes

- 1. Prepare the following stock solutions:
 - 200 mM THPTA ligand in water
 - 100 mM CuSO4 in water
 - Azide-modified peptide in water or DMF (depending on your peptide solubility, >10 mg/ml if possible)
 - 100 mM sodium ascorbate in water
 - 10 mM dye alkyne in DMSO or water (see our website for recommended solvent)
- 2. Incubate CuSO₄ with THPTA ligand in a 1:2 ratio several minutes before the reaction. This solution is stable for several weeks when frozen.
- 3. To the azide-modified peptide solution, add an excess of dye alkyne (5-10equivalents by molar ratio).
- 4. Add 5-10 equivalents of THPTA/CuSO₄.
- 5. Add 10-20 equivalents of sodium ascorbate.
- 6. Stir, vortex or shake the reaction mixture at room temperature for 30-60 minutes.
- 7. Purify your desired peptide by HPLC.

Labeling Small Organic Alkyne Molecules with Dye Alkynes

- 1. Prepare the following stock solutions:
 - 200 mM THPTA ligand in water
 - 100 mM CuSO4 in water
 - Alkyne compound in water or DMF (depending on your compound solubility, >10 mg/ml if possible,)
 - 100 mM sodium ascorbate in water
 - 10 mM dye alkyne in DMSO or water (see our website for recommended solvent).
- 2. Incubate CuSO₄ with THPTA ligand in a 1:2 ratio several minutes before the reaction. This solution is stable for several weeks when frozen.
- 3. To the alkyne solution, add an excess of dye alkyne (5-10 equivalents by molar ratio).
- 4. Add 25 equivalents of THPTA/CuSO₄.
- 5. Add 50 equivalents of sodium ascorbate.
- 6. Stir the reaction mixture at room temperature for 30-60 minutes.
- 7. Purify your desired molecule by chromatography or other methods.

Labeling Biopolymers with Dye Alkynes

- 1. Prepare the following stock solutions:
 - 200 mM THPTA ligand in water
 - 100 mM CuSO4 in water
 - Azide-modified biopolymer in water (as concentrated as possible, e.g., >5 mg/ml)
 - 100 mM sodium ascorbate in water
 - 10 mM dye alkyne in DMSO or water (see our website for recommended solvent).
- 2. Incubate CuSO₄ with THPTA ligand in a 1:2 ratio several minutes before the reaction. This solution is stable for several weeks when frozen.
- 3. To the azide-modified biopolymer solution, add an excess of dye alkyne (Loading ratio: 5-20 dye alkyne/alkyne).
- 4. Add 5 molar equivalents (referenced to dye alkyne) of THPTA/CuSO₄.
- 5. Add 10 equivalents of sodium ascorbate (referenced to dye alkyne).
- 6. Stir, vortex or shake the reaction mixture at room temperature for 30-60 minutes.
- 7. Purify your desired molecule by gel filtration or dialysis.

Labeling Cells, Cell Lysates or Biological Samples with Dye alkynes or Dye Alkynes

- 1. Prepare the following click solutions:
 - 100 mM THPTA ligand in aqueous buffer or water
 - 20 mM CuSO4 in water
 - 300 mM sodium ascorbate in water
 - 2.5 mM alkyne or azide labeling reagent in water or DMSO
- 2. For each azide- or azide-modified cell or cell lysate sample, add the following reagents to a 1.5 mL microfuge tube, then vortex briefly to mix.
 - 50 µL cell or cell lysate sample
 - 50 µL PBS buffer
 - 50 μL of 5 mM corresponding dye alkyne (or dye alkyne) detection reagent in DMSO or water
- 3. Add $10 \,\mu\text{L}$ of $100 \,\text{mM}$ THPTA solution, vortex briefly to mix.
- 4. Add 10 μL of 20 mM CuSO4 solution, vortex briefly to mix.
- 5. Add 10 μ L of 300 mM sodium ascorbate solution to initiate click reaction, vortex briefly to mix.
- 6. Protect reaction from light and allow click reaction to incubate for 30 minutes at room temperature.
- 7. Cells or cell lysates are now click labeled and ready for downstream processing and/or analysis.