

## Blunting DNA ends by 3'-overhang removal or fill-in of 3' recessed-end Utilizing T4 DNA Polymerase

### Introduction

Polymerase Chain Reaction (PCR) is a powerful technique used to amplify DNA through the use of the enzyme T4 DNA Polymerase. GoldBio T4 DNA Polymerase has DNA-dependent DNA polymerase activity and potent 3'→5' exonuclease activity. GoldBio T4 DNA Polymerase is ideal for 3'-overhang removal or 5'-overhang fill-in to form blunt ends, probe labeling using replacement synthesis, DNA library preparation for Next-generation sequencing, ligation-independent cloning of PCR products and second strand synthesis in site-directed mutagenesis. Here, we describe a general protocol for the use of GoldBio T4 DNA Polymerase in the blunting of ends by 3' overhang removal or 3' recessed-end fill-in.

### Materials

- T4 DNA Polymerase (GoldBio Catalog # [T-412](#))
- 10x T4 DNA Polymerase Buffer

Not supplied:

- 100µM dNTPs (GoldBio Catalog # [D-900](#))
- H<sub>2</sub>O PCR Grade

**Note:** The storage buffer contains 50mM Tris-HCl, 50mM KCl, 1mM DTT, 0.1mM EDTA, 50% glycerol, pH 7.5 at 25°C.

**Note:** 10x T4 DNA Polymerase Buffer contains 500mM Tris-HCl, 100mM MgCl<sub>2</sub>, 50mM DTT, pH 7.5 at 25°C.

**Note:** T4 DNA Polymerase concentration is 5 units/ul. One unit of T4 DNA Polymerase converts 10 nmol of dNTPs into acid-insoluble material in 30 minutes at 37°C under standard assay conditions.

**Note:** T4 DNA Polymerase is inactivated by incubating at 70°C for 15 minutes.

### Storage and Handling

- Store GoldBio T4 DNA Polymerase and the 10x T4 DNA Polymerase Buffer at -20°C.
- These products may be shipped on blue ice and should be stored at -20°C immediately upon arrival. When stored under the recommended conditions and handled correctly, these products should be stable for at least 1 year from the date of receipt.

- Thaw on ice and mix by gentle vortexing. After thawing, these products should be kept on ice before use. These products can be refrozen for storage.

## Method

Blunting ends by 3' overhang removal or 3' recessed-end fill-in.

1. Dissolve linear DNA or PCR product in 1x reaction buffer supplemented with 100 $\mu$ M dNTPs.
2. Add 1 unit T4 DNA Polymerase per  $\mu$ g DNA.
3. Incubate at room temperature for 5-30 minutes.
4. Stop the reaction by heating at 70°C for 20 minutes.

## Associated Products

- [dNTP Mix \(GoldBio Catalog # D-900\)](#)
- [T4 Polynucleotide Kinase \(PNK\) \(GoldBio Catalog # T-413\)](#)
- [T4 UvsX DNA Recombinase \(GoldBio Catalog # T-414\)](#)
- [T4 UvsY Protein \(GoldBio Catalog # T-415\)](#)
- [T4 gp32 Protein \(GoldBio Catalog # T-416\)](#)
- [T4 DNA Helicase \(gp41\) \(GoldBio Catalog # T-417\)](#)
- [T4 gp46 Protein \(GoldBio Catalog # T-418\)](#)
- [T4 gp47 Protein \(GoldBio Catalog # T-419\)](#)
- [T4 gp59 Protein \(GoldBio Catalog # T-420\)](#)
- [T4 DNA Ligase – 400 units/ \$\mu\$ l \(GoldBio Catalog # T-410\)](#)
- [T4 DNA Ligase – 2000 units/ \$\mu\$ l \(GoldBio Catalog # T-411\)](#)

## References

- Tabor, S. and Struhl, K. 1989. DNA-dependent DNA polymerases. In F. M. Ausebel, R. Brent, R. E. Kingston, D. D. Moore, J. G. Seidman, J. A. Smith and K. Struhl (Ed.). *Current Protocols in Molecular Biology*, 3.5.10-3.5.12. New York: John Wiley & Sons, Inc.
- Sambrook, J., Fritsch, E. F. and Maniatis, T. 1989. *Molecular Cloning: A Laboratory Manual*. (2<sup>nd</sup> ed), 5.44-5.47. Cold Spring Harbor: Cold Spring Harbor Laboratory Press.
- Kunkel, T. A., Roberts, J. D. and Zakour, R. A. 1987. Rapid and efficient site-specific mutagenesis without phenotypic selection. *Methods Enzymol*, 154:367-82.