

Ligation Protocol Utilizing T4 DNA Ligase

Introduction

GoldBio T4 DNA Ligase catalyzes the formation of a phosphodiester bond between juxtaposed 5'-phosphate and 3'-hydroxyl termini in duplex DNA or RNA. GoldBio T4 DNA ligase joins DNA fragments with either cohesive or blunt termini and repairs single-stranded nicks in duplex DNA, RNA or DNA/RNA hybrids. In addition, GoldBio T4 DNA ligase can be used to clone restriction enzyme-generated DNA fragments and PCR products. T4 DNA ligase can also be used for Next-gen library preparation and self-circularization of linear DNA. Finally, T4 DNA ligase can also join linkers and adapters to cohesive or blunt-ended DNA and repair nicks in duplex DNA, RNA, or DNA/RNA hybrids. Here, we describe a general protocol for the use of GoldBio T4 DNA Ligase to ligate vector and insert DNA.

Materials

- T4 DNA Ligase (GoldBio Catalog # [T-410](#) or [T-411](#))
- 10x T4 DNA Ligase Reaction Buffer

Not supplied:

- H₂O PCR grade

Note: T4 DNA Ligase is inhibited by metal chelators, phosphate and ammonium ions, KCl and NaCl at a concentration higher than 50mM. It can be inactivated by incubating at 70°C for 15 minutes or by adding EDTA.

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

Note: 1x T4 DNA Ligase Reaction Buffer contains 50mM Tris-HCl, 10mM MgCl₂, 10mM DTT, 1mM ATP, pH 7.5 at 25°C.

Note: One unit is defined as the amount of enzyme required to give 50% ligation of HindIII fragments of λ DNA (250 ng/ μ l) in a total reaction volume of 20 μ l in 30 minutes at 16°C in 1x T4 DNA ligase reaction buffer.

Storage and Handling

- Store GoldBio T4 DNA Ligase and the 10x T4 DNA Ligase Reaction 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

1. Set up a reaction according to Table 1, in a microcentrifuge tube, on ice. Use a molar ratio of 1:3 vector to insert DNA (see Calculations section).

Table 1. Set up of ligation reaction (400 units/ μ l Ligase)

Component	20 μ l Reaction
Vector DNA	x μ l
Insert DNA	x μ l
10x T4 Ligase Buffer	2.0 μ l
T4 DNA Ligase	1.0 μ l
Nuclease-free water	Up to 20.0 μ l

2. Gently mix the reaction and centrifuge briefly.
3. For cohesive ends, incubate at room temperature for 10 minutes or 16°C overnight.
4. For blunt ends, incubate at room temperature for 2 hours or 16°C overnight.
5. Heat-inactivate at 70°C for 15 minutes.
6. Cool on ice and transform 2 μ l of the reaction into 50 μ l of competent cells.

Calculations

1. To calculate ng of insert to add to ligation reaction with known amount of vector, first calculate the ratio coefficient (x).

$$\frac{bp \text{ size vector}}{bp \text{ size insert}} = x$$

2. Determine the quantity of vector (y) to use.

Note: Generally, 10 ng of vector should be sufficient.

3. Calculate the quantity of insert to use at 1:1 ratio (z).

$$\frac{y}{x} = z$$

4. Calculate the quantity of insert needed for a 3:1 insert:vector molar ratio ligation reaction.

$$3 \times z = \text{Quantity of insert}$$

Associated Products

- [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 – 2000 units/μl \(GoldBio Catalog # T-411\)](#)

References

Engler, M. J. and Richardson, C. C. 1982. DNA Ligases. In: P.D. Boyer (Ed.) The Enzymes, vol. XV, Academic Publishers, New York, pp. 3-29.