



**T4 DNA Ligase**

*E. coli* lambda lysogen NM 989

| Cat. No. | Amount                                   |
|----------|--|
| EN-149S  | 400 Weiss units (80000 CE units)         |
| EN-149L  | 5 x 400 Weiss units (5 x 80000 CE units) |

**Unit Definition:** One Weiss unit is defined as the amount of enzyme required to catalyze the exchange of 1 nmol of <sup>32</sup>P from pyrophosphate to ATP, into Norit-adsorbable material in 20 minutes at 37 °C.

**For general laboratory use.**

**Shipping:** shipped on gel packs

**Storage Conditions:** store at -20 °C

**Additional Storage Conditions:** avoid freeze/thaw cycles

**Shelf Life:** 12 months

**Form:** liquid (Supplied in 10 mM Tris-HCl pH 7.4, 50 mM KCl, 0.1 mM EDTA, 1 mM DTT, 200 µg/ml BSA and 50 % [v/v] glycerol)

**Concentration:** 2.5 Weiss units/µl (500 CE units/µl)

**Description:**

T4 DNA Ligase catalyzes the formation of a phosphodiester bond between juxtaposed 5' phosphate and 3'-hydroxyl termini in duplex DNA or RNA.

**Content:**

**Standard Ligation Buffer, 10x conc.**

500 mM Tris-HCl pH 7.8 at 25 °C, 100 mM MgCl<sub>2</sub>, 100 mM DTT, 10 mM ATP and 25 µg/ml BSA

**Fast Ligation Buffer, 2x conc.**

60 mM Tris-HCl pH 7.8 at 25 °C, 20 mM MgCl<sub>2</sub>, 20 mM DTT, 2 mM ATP and 10 % PEG

| component                           | EN-149S | EN-149L    |
|-------------------------------------|---------|------------|
| T4 DNA Ligase                       | 160 µl  | 5 x 160 µl |
| Standard Ligation Buffer, 10x conc. | 1 ml    | 5 x 1 ml   |
| Fast Ligation Buffer, 2x conc.      | 5 ml    | 5 x 5 ml   |

**Heat inactivation:**

T4 DNA Ligase can be inactivated by incubation at 65 °C for 10 minutes.

**Note:**

- One Cohesive-End Ligation Unit (CEU) is defined as the amount of enzyme required to give 50 % ligation of *Hind* III fragments of λ DNA (5' DNA termini concentration of 0.12 µM, 300 µg/ml) in a total reaction volume of 20 µl in 30 minutes at 16 °C in 1x T4 DNA Ligase Reaction Buffer.
- One Weiss unit is equivalent to approx. 200 CE units.
- T4 DNA Ligase is strongly inhibited by NaCl or KCl if the concentration exceeds 200 mM.
- Ligation of blunt-ended and single-base pair overhang fragments requires about 50 times as much enzyme to achieve the same extent of ligation as cohesive-end DNA fragments. Blunt-end ligation may be enhanced by addition of PEG 4000 (10 % w/v final concentration) or hexamine chloride, or by reducing the ATP concentration to 50 µM.
- To dilute T4 DNA Ligase for subsequent storage at -20 °C a storage buffer containing 50 % glycerol should be used, to dilute Ligase for immediate use, 1x Reaction Buffer is recommended.



## T4 DNA Ligase

*E. coli* lambda lysogen NM 989

### Assay Set-Up:

#### Standard Ligation Assay

| comp.                               | final amount/conc.  | 20 µl assay      |
|-------------------------------------|---------------------|------------------|
| Standard Ligation Buffer, 10x conc. | 1x                  | 2 µl             |
| Vector/Insert DNA                   | 100 ng - 1 µg       | 100 ng - 1 µg    |
| T4 DNA Ligase                       | 0.1 - 1 Weiss units | 0.04-0.4 µl      |
| PCR-grade Water                     | -                   | fill up to 20 µl |

Incubate for 20 - 30 min at 16 °C for optimal ligation.

#### Fast Ligation Assay

| comp.                          | final amount/conc.  | 20 µl assay      |
|--------------------------------|---------------------|------------------|
| Fast Ligation Buffer, 2x conc. | 1x                  | 10 µl            |
| Vector/Insert DNA              | 100 ng - 1 µg       | 100 ng - 1 µg    |
| T4 DNA Ligase                  | 0.1 - 1 Weiss units | 0.04-0.4 µl      |
| PCR-grade Water                | -                   | fill up to 20 µl |

Incubate for 5 min for cohesive-ended ligations or 15 min for blunt-ended ligations at ambient temperature (20 - 25 °C).