

## T4 DNA Ligase (Cloned)

Store at  $-20^{\circ}\text{C}$ .  
*Do not store in a frost-free freezer.*

|                                       |  |
|---------------------------------------|--|
| <b>Catalog #:</b>                     | AM2134   |
| <b>Product Description:</b>           | High-purity T4 DNA Ligase, for use in ligation of blunt- and cohesive-ended DNA, and for nick repair in duplex DNA.                                    |
| <b>Source:</b>                        | An <i>E. coli</i> strain overexpressing T4 DNA Ligase.   |
| <b>Amount:</b>                        | 1000 Units   |
| <b>Unit Concentration:</b>            | 5 U/ $\mu\text{L}$ (Weiss)   |
| <b>Unit Definition:</b>               | One unit catalyzes the exchange of 1 nmol of inorganic phosphate from pyrophosphate into Norit-adsorbable material in 20 min at $37^{\circ}\text{C}$ . |
| <b>Additional Materials Included:</b> | 500 $\mu\text{L}$ 10X T4 DNA Ligase Reaction Buffer<br>500 mM Tris, pH 7.8<br>100 mM $\text{MgCl}_2$<br>100 mM DTT<br>10 mM ATP<br>0.25 mg/mL BSA      |

*Note: Precipitates may form during storage; warm to  $37^{\circ}\text{C}$  and vortex to resuspend any precipitated material.*

|                            |   |
|----------------------------|---|
| <b>Storage Conditions:</b> | Store at $-20^{\circ}\text{C}$ . <b>Do not store in a frost-free freezer.</b>               |
| <b>Storage Buffer:</b>     | (Not included) 20 mM Tris-HCl (pH 7.6), 60 mM KCl, 1 mM DTT, 1 mM EDTA, 50% glycerol (v/v). |

---

### USER INFORMATION

---

|                             |  |
|-----------------------------|--|
| <b>General Information:</b> | When incubated in the presence of ATP (included in the supplied 10X Reaction Buffer), T4 DNA Ligase (E.C. 6.5.1.1) catalyzes the formation of phosphodiester bonds between the 5'-phosphoryl group and the 3'-hydroxyl group of double-stranded DNA fragments. T4 DNA Ligase will also close single-stranded "nicks" in double-stranded DNA molecules.   |
| <b>Applications:</b>        | The following general protocol is adapted from <i>Current Protocols in Molecular Biology</i> .<br><br>Mix together:<br>5 $\mu\text{L}$ 10X T4 DNA Ligase Reaction Buffer (included)<br>1 $\mu\text{g}$ DNA<br>1–5 U T4 DNA Ligase<br>To 50 $\mu\text{L}$ with nuclease-free water<br><br>For cohesive (sticky) end ligations, incubate at room temperature for 1–3 hr. If poor results are obtained with a 1–3 hr incubation, incubating overnight at $16^{\circ}\text{C}$ may improve the ligation reaction.<br><br>For blunt end ligation, incubate overnight at $16^{\circ}\text{C}$ .<br><br>The reaction mixture can be used for transformation without further processing. |
| <b>Reference:</b>           | Ausubel FM, Brent R, Kingston RE, Moore DD, Seidman JG, Smith JA, Struhl K, editors. (2006) Enzymatic Manipulation of DNA and RNA (Chapter 3). In <i>Current Protocols in Molecular Biology</i> . John Wiley & Sons, Inc.  |

---

### QUALITY CONTROL

---

T4 DNA Ligase is tested for contaminating nonspecific endonuclease, exonuclease, and protease activity. Functionally is determined in a ligation reaction with HindIII-cut Lambda DNA.

---

### OTHER INFORMATION

---

|                                     |   |
|-------------------------------------|---|
| <b>Material Safety Data Sheets:</b> | Material Safety Data Sheets (MSDSs) can be printed or downloaded from product-specific links on our website at the following address: <a href="http://www.ambion.com/techlib/msds">www.ambion.com/techlib/msds</a> . Alternatively, e-mail your request to <a href="mailto:MSDS_Inquiry_CCRM@appliedbiosystems.com">MSDS_Inquiry_CCRM@appliedbiosystems.com</a> . Specify the catalog or part number(s) of the product(s), and we will e-mail the associated MSDSs unless you specify a preference for fax delivery. For customers without access to the internet or fax, our technical service department can fulfill MSDS requests placed by telephone or postal mail. (Requests for postal delivery require 1–2 weeks for processing.) |
|-------------------------------------|---|

**Warranty and Liability:**

*For research use only. Not for use in diagnostic procedures.*

Applied Biosystems is committed to delivering superior product quality and performance, supported by industry-leading global service and technical support teams. Warranty information for the accompanying consumable product is available at [www.ambion.com/info/warranty](http://www.ambion.com/info/warranty) in "Limited Warranty for Consumables," which is subject to the exclusions, conditions, exceptions, and limitations set forth under the caption "EXCLUSIONS, CONDITIONS, EXCEPTIONS, AND LIMITATIONS" in the full warranty statement. Please contact Applied Biosystems if you have any questions about our warranties or would like information about post-warranty support.

**Trademarks, Patents, and Licensing:**

Applied Biosystems, AB (Design), and Ambion are registered trademarks of Applera Corporation or its subsidiaries in the US and/or certain other countries. All other trademarks are the sole property of their respective owners.

© 2007 Ambion, Inc. All rights reserved. 4393895A