

## Product Contents

### Shrimp Alkaline Phosphatase (SAP), Blue/White Cloning Qualified:

| Part No. | Conc. (u/μl) | Size (units) |
|----------|--------------|--------------|
| M820A    | 1            | 500          |

**Description:** Shrimp Alkaline Phosphatase (SAP) catalyzes the dephosphorylation of 5' phosphates from DNA. Unlike Calf Intestinal Alkaline Phosphatase, SAP is completely and irreversibly inactivated by heating at 65°C for 15 minutes. Alkaline phosphatases are used to prevent recircularization and religation of linearized cloning vehicle DNA by removing phosphate groups from both 5'-termini and may also be used for the dephosphorylation of 5' phosphorylated ends of DNA for subsequent labeling with [<sup>32</sup>P]ATP and T4 Polynucleotide Kinase. SAP is active on 5' overhangs, 5' recessed and blunt ends (1).

**Enzyme Storage Buffer:** Shrimp Alkaline Phosphatase (SAP) is supplied in 25mM Tris-HCl (pH 7.6 at 4°C), 1mM MgCl<sub>2</sub>, 0.1mM ZnCl<sub>2</sub> and 50% glycerol.

**Shrimp Alkaline Phosphatase 10X Reaction Buffer (M821A):** When the 10X Reaction Buffer supplied with this enzyme is diluted 1:10, it has a composition of 50mM Tris-HCl (pH 9.0 at 37°C) and 10mM MgCl<sub>2</sub>.

**Source:** Shrimp (*Pandalus borealis*).

**Unit Definition:** One unit is defined as the amount of enzyme required to catalyze the hydrolysis of 1μmol of 4-nitrophenyl phosphate per minute at 37°C in 1M diethanolamine, 10.9mM 4-nitrophenyl phosphate, 0.5mM MgCl<sub>2</sub> (pH 9.8). See the unit concentration on the Product Information Label.

**Storage Temperature:** See the Product Information Label for storage recommendations. Avoid multiple freeze-thaw cycles. See the expiration date on the Product Information Label.

Part# 9PIM820

Revised 3/07



AF9P1 M820 0307M820



**Promega**

#### Promega Corporation

|                        |  |
|------------------------|--|
| 2800 Woods Hollow Road |  |
| Madison, WI 53711-5399 | USA  |
| Telephone              | 608-274-4330   |
| Toll Free              | 800-356-9526   |
| Fax                    | 608-277-2516   |
| Internet               | <a href="http://www.promega.com">www.promega.com</a> |

## Quality Control Assays

### Contaminant Activity

**Endonuclease Assay:** To test for endonuclease activity, 1μg of Type I supercoiled plasmid DNA is incubated with 5 units of Shrimp Alkaline Phosphatase in 1X Reaction Buffer for one hour at 37°C. Following incubation, the supercoiled DNA is visualized on an ethidium bromide-stained agarose gel to verify the absence of visible nicking or cutting.

**DNase and RNase Assay:** To test for nuclease activity, 50ng of radiolabeled DNA or radiolabeled RNA is incubated with 5 units of Shrimp Alkaline Phosphatase in 1X Reaction Buffer for one hour at 37°C, and the release of radiolabeled nucleotides is monitored by scintillation counting of TCA-soluble material. Minimum passing specification is ≤3% release for DNase and ≤3% release for RNase.

**Blue/White Assay:** pGEM®-3Zf(+) Vector is digested with representative restriction enzymes (leaving 5'-termini, 3'-termini or blunt ends). Each microgram of cut plasmid is treated with 2 units of Shrimp Alkaline Phosphatase for 1 hour at 37°C, kinased and ligated. The religated plasmid is then transformed into JM109 cells that are plated on X-Gal/IPTG/Ampicillin plates. White colonies result from transformation with ligated plasmids with damaged ends. These white colonies represent the number of false positives expected in a typical cloning experiment. Enzymes that generate overhangs must produce fewer than 2% white colonies, and blunt-cutting enzymes must produce fewer than 5% white colonies (2).

## References

1. Sambrook, J., Fritsch, E.F. and Maniatis, T. (1989) *Molecular Cloning: A Laboratory Manual*, Cold Spring Harbor Laboratory Press, Cold Spring Harbor, New York.
2. Hung, L. *et al.* (1991) A blue/white cloning assay for quality control of DNA restriction and modifying enzymes. *Promega Notes* **33**, 12-13.

#### PRODUCT USE LIMITATIONS, WARRANTY, DISCLAIMER

Promega manufactures products for a number of intended uses. Please refer to the product label for the intended use statements for specific products. Promega products contain chemicals which may be harmful if misused. Due care should be exercised with all Promega products to prevent direct human contact.

Each Promega product is shipped with documentation stating specifications and other technical information. Promega products are warranted to meet or exceed the stated specifications. Promega's sole obligation and the customer's sole remedy is limited to replacement of products free of charge in the event products fail to perform as warranted. Promega makes no other warranty of any kind whatsoever, and SPECIFICALLY DISCLAIMS AND EXCLUDES ALL OTHER WARRANTIES OF ANY KIND OR NATURE WHATSOEVER, DIRECTLY OR INDIRECTLY, EXPRESS OR IMPLIED, INCLUDING, WITHOUT LIMITATION, AS TO THE SUITABILITY, PRODUCTIVITY, DURABILITY, FITNESS FOR A PARTICULAR PURPOSE OR USE, MERCHANTABILITY, CONDITION, OR ANY OTHER MATTER WITH RESPECT TO PROMEGA PRODUCTS. In no event shall Promega be liable for claims for any other damages, whether direct, incidental, foreseeable, consequential, or special (including but not limited to loss of use, revenue or profit), whether based upon warranty, contract, tort (including negligence) or strict liability arising in connection with the sale or the failure of Promega products to perform in accordance with the stated specifications.

© 2000-2007 Promega Corporation. All Rights Reserved.

pGEM and Wizard are registered trademarks of Promega Corporation. MULTICORE is a trademark of Promega Corporation.

Products may be covered by pending or issued patents or may have certain limitations. Please visit our Web site for more information.

All specifications are subject to change without prior notice.

Product claims are subject to change. Please contact Promega Technical Services or access the Promega online catalog for the most up-to-date information on Promega products.

Part# 9PIM820

Printed in USA. Revised 3/07

## I. Reaction Conditions

### A. Standard Dephosphorylation and Ligation Reaction

1. Incubate SAP (1 unit/ $\mu$ g DNA) with restriction-digested vector (restriction enzyme and buffer removed, DNA in water or TE buffer) at 37°C for 15 minutes in 1X SAP reaction buffer in a final volume of 30–50 $\mu$ l. This is a sufficient amount of SAP to completely dephosphorylate the vector regardless of overhang type (5', 3' or blunt).
2. Inactivate SAP by heating to 65°C for 15 minutes.
3. Centrifuge the reaction mix. Remove a 1–2 $\mu$ l aliquot for ligation with appropriate DNA insert in 1X ligation buffer at 4°C overnight for all types of DNA overhangs. Final volume should be between 10–50 $\mu$ l. Molar ratios for vector to insert can range between 1:8 to 8:1. A 1:1 ratio is recommended as a starting point.
4. Transform the ligated material directly into competent *E. coli* cells.

### B. Streamlined Restriction Digestion, Dephosphorylation and Ligation Procedure

1. Combine restriction digestion and dephosphorylation of DNA vector in 1X restriction enzyme buffer. Use 15 units of restriction enzyme/ $\mu$ g vector and 10 units SAP/ $\mu$ g vector in a final volume of 30–50 $\mu$ l. Incubate at 37°C for 15 minutes. This is a sufficient amount of SAP to completely dephosphorylate the vector regardless of overhang type (5', 3' or blunt).
2. Heat-inactivate both restriction enzyme and SAP for 15 minutes at 65°C.  
**Note:** Not all restriction enzymes can be heat inactivated.
3. Centrifuge and remove 1–2 $\mu$ l (~40ng) of vector for ligation with appropriate DNA insert using 2X Rapid Ligation Buffer (Cat.# C6711) and 2 $\mu$ l (6 units) of T4 DNA Ligase (Cat.# M1801) at 15°C for 5 minutes (3' or 5' ends) or 15 minutes for blunt ends in a final reaction volume of 10–50 $\mu$ l. We recommend starting with a 1:2 molar ratio of vector:insert DNA.
4. Transform the ligated material directly into competent *E. coli* cells.

### C. Standard Dephosphorylation and Labeling Reaction Using T4 Polynucleotide Kinase

1. Incubate SAP (1 unit/0.5pmol ends) and DNA at 37°C for 15 minutes in 1X SAP buffer in a final volume of 30–100 $\mu$ l.
2. Purify the DNA using the Wizard® DNA Clean-Up System (Cat.# A7280).
3. Labeling is carried out by incubating 8–10 units T4 Polynucleotide Kinase (M4101),  $\leq$ 10pmol DNA ends and 15 $\mu$ l [ $\gamma$ -<sup>32</sup>P] ATP (3,000Ci/mmol, 10mCi/ml) at 37°C for 10 minutes in 1X PNK buffer and a final volume of 50 $\mu$ l.

## II. Composition of Buffers and Solutions

### 2X Rapid Ligation Buffer

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

## III. Related Products

| Product                     | Size            | Cat.# |
|-----------------------------|-----------------|-------|
| Wizard® DNA Clean-Up System | 100 preps       | A7280 |
| T4 Polynucleotide Kinase    | 100u            | M4101 |
| 2X Rapid Ligation Buffer    | 3 × 500 $\mu$ l | C6711 |
| T4 DNA Ligase               | 1,000u          | M1801 |

**Table 1. Activity of Shrimp Alkaline Phosphatase (Part# M820A) in Restriction Enzyme Buffers vs. 1X Shrimp Alkaline Phosphatase Buffer.**

| Promega Restriction Enzyme Buffer | % Activity of SAP |
|-----------------------------------|-------------------|
| A                                 | 20                |
| B                                 | 20                |
| C                                 | 25                |
| D                                 | 35                |
| E                                 | 20                |
| F                                 | 60                |
| G                                 | 30                |
| H                                 | 30                |
| J                                 | 25                |
| K                                 | 20                |
| L                                 | 30                |
| MULTI-CORE™                       | 10                |