

## The Signature Series

### Streptavidin Alkaline Phosphatase Detection System

Catalog no. 750144, 25 ml

---

**Reagent Design**

The Streptavidin Alkaline Phosphatase (AP) Detection System combines the stability and high sensitivity of alkaline phosphatase with the specificity of streptavidin. The Streptavidin AP Detection System combines with the biotin moieties in the Universal Secondary Antibody or in the probe. This combination localizes the *in situ* immune interaction of the primary antibody and the tissue antigen or of the probe and the probe target. The reagent contains optimally diluted, high specific activity streptavidin conjugated alkaline phosphatase diluted in the Detection System Diluent and is stable for six months when stored at 4 °C. This reagent is color coded orange for easy identification.

---

**Reagent Use**

In immunoalkaline phosphatase and *in situ* hybridization procedures with Stable Naphthol Phosphate and Stable Fast Red binary chromogen system, the Streptavidin AP Detection System generates a permanent, red immunostain. Applying the Streptavidin AP Detection System at 33-50 °C for four minutes yields sensitive and uniform immunostaining. Protocols for using this reagent in MicroProbe™ immunocytochemistry and *in situ* hybridization are available from Invitrogen.

---

**Quality Control**

This detection system has been thoroughly tested and verified for immunocytochemistry and *in situ* hybridization using standard, manual technology and a MicroProbe™ capillary action system. In addition, it has been thoroughly tested and verified for immunocytochemistry using a CodeOn™ Automated Molecular Pathology System.

---

**Signature**

The Streptavidin AP Detection System was designed and developed by David J. Brigati, M. D. with the assistance of our staff.

---

**Technical Service**

For more information or technical assistance, call, write, fax, or email. Additional international offices are listed on our Web page ([www.invitrogen.com](http://www.invitrogen.com)).

**Corporate Headquarters:**

Invitrogen Corporation  
1600 Faraday Avenue  
Carlsbad, CA 92008 USA  
Tel: 1 760 603 7200  
Tel (Toll Free): 1 800 955 6288  
Fax: 1 760 602 6500  
E-mail:  
[tech\\_service@invitrogen.com](mailto:tech_service@invitrogen.com)

**Japanese Headquarters:**

Invitrogen Japan K.K.  
Nihonbashi Hama-Cho Park Bldg. 4F  
2-35-4, Hama-Cho, Nihonbashi  
Tel: 81 3 3663 7972  
Fax: 81 3 3663 8242  
E-mail: [jpinfo@invitrogen.com](mailto:jpinfo@invitrogen.com)

**European Headquarters:**

Invitrogen Ltd  
Inchinnan Business Park  
3 Fountain Drive  
Paisley PA4 9RF, UK  
Tel: +44 (0) 141 814 6100  
Tel (Toll Free): 0800 5345 5345  
Fax: +44 (0) 141 814 6117  
E-mail: [eurotech@invitrogen.com](mailto:eurotech@invitrogen.com)

---

©2001-2002 Invitrogen Corporation. All rights reserved.

MicroProbe™ is a trademark of the Fisher Scientific Co., Pittsburgh, PA. CodeOn™ is a trademark of Instrumentation Laboratory, Lexington, MA.