

The Signature Series

Stable Fast Red TR/Stable Naphthol Phosphate

Catalog no. 750152; 125 ml of each

Reagent Design

Stable Fast Red TR / Stable Naphthol Phosphate is an easy to use, stable, sensitive, low cost, high volume, prediluted chromogen system for immunoalkaline phosphatase and *in situ* hybridization procedures. Its localization is unambiguous and without background, and it creates a permanent record of the results under Pristine Mount without fading. Stable Fast Red TR and Stable Naphthol Phosphate work together to produce a sensitive, intense, red chromogen for alkaline phosphatase in both immunocytochemistry and *in situ* hybridization. Both reagents are stable when stored separately at 4 °C for one year.

Reagent Use

Stable Fast Red TR and Stable Naphthol Phosphate are used together in immunoalkaline phosphatase and *in situ* DNA hybridization procedures to localize genes, their transcripts, and protein antigen products. To prepare the working chromogen, pour any volume of Stable Fast Red TR into a graduated conical centrifuge tube, and add an equal volume of Stable Naphthol Phosphate. For optimal signal enhancement, use the working chromogen immediately after mixing the Stable Fast Red TR and the Stable Naphthol Phosphate solutions together. Only two 10-minute, 45 °C applications of the combined working Stable Fast Red TR / Stable Naphthol Phosphate chromogen are necessary to produce powerful, permanent, and clean results. However, up to four additional applications of this binary chromogen can increase immunostaining. Mount tissue sections in Pristine Mount for permanent preservation. Do not dehydrate tissues with alcohol or xylene after staining or use a xylene-based mounting media, because xylene and alcohol eliminate the chromogen signal. Instead, dry the tissues at room temperature after the final staining step, and then mount them with a water-based mounting medium, such as Pristine Mount, either with or without a cover slip.

Quality Control

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

Signature

Stable Fast Red TR / Stable Naphthol Phosphate was designed and developed by David J. Brigati, M. D. with the assistance of our staff.

Continued on next page

Stable Fast Red TR/Stable Naphthol Phosphate, Continued

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).

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

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

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

©2001-2002 Invitrogen Corporation. All rights reserved.

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