



00-2209 1 Kit
BCIP/NBT Kit
Lot No.

BCIP/NBT SUBSTRATE KIT

SUBSTRATE FOR ALKALINE PHOSPHATASE

FORM: liquid.

KIT CONTENTS:

- A - Concentrated BCIP/Buffer Solution (10x), 100 ml
- B - Concentrated NBT solution (10x), 100 ml

BCIP - 5-Bromo-4-Chloro-3-Indolyl phosphate
NBT - Nitroblue tetrazolium salt

BLOTTING PROCEDURE:

BCIP is an insoluble substrate for alkaline phosphatase which is especially good for staining blots.

Add 1 ml of each reagent to 8 ml of distilled water immediately before use. Blots will normally require 10-20 ml. Incubate the blots in this solution at room temperature for 10-30 min. until color develops. Do not overdevelop as this may produce background. Discard the unused portion of the diluted substrate solution.

STORAGE:

Store at 2-8°C.

REMARKS:

For research use only. Not for human or drug use. No preservative added.

www.invitrogen.com

Invitrogen Corporation • 542 Flynn Rd • Camarillo • CA 93012 • Tel: 800.955.6288
E-mail: techsupport@invitrogen.com

PI002209

(Rev 10/08) DCC-08-1089

Important Licensing Information - These products may be covered by one or more Limited Use Label Licenses (see the Invitrogen Catalog or our website, www.invitrogen.com). By use of these products you accept the terms and conditions of all applicable Limited Use Label Licenses. Unless otherwise indicated, these products are for research use only and are not intended for human or animal diagnostic, therapeutic or commercial use.



00-2209 1 Kit
BCIP/NBT Kit
Lot No.

BCIP/NBT SUBSTRATE KIT

SUBSTRATE FOR ALKALINE PHOSPHATASE

FORM: liquid.

KIT CONTENTS:

- A - Concentrated BCIP/Buffer Solution (10x), 100 ml
- B - Concentrated NBT solution (10x), 100 ml

BCIP - 5-Bromo-4-Chloro-3-Indolyl phosphate
NBT - Nitroblue tetrazolium salt

BLOTTING PROCEDURE:

BCIP is an insoluble substrate for alkaline phosphatase which is especially good for staining blots.

Add 1 ml of each reagent to 8 ml of distilled water immediately before use. Blots will normally require 10-20 ml. Incubate the blots in this solution at room temperature for 10-30 min. until color develops. Do not overdevelop as this may produce background. Discard the unused portion of the diluted substrate solution.

STORAGE:

Store at 2-8°C.

REMARKS:

For research use only. Not for human or drug use. No preservative added.

www.invitrogen.com

Invitrogen Corporation • 542 Flynn Rd • Camarillo • CA 93012 • Tel: 800.955.6288
E-mail: techsupport@invitrogen.com

PI002209

(Rev 10/08) DCC-08-1089

Important Licensing Information - These products may be covered by one or more Limited Use Label Licenses (see the Invitrogen Catalog or our website, www.invitrogen.com). By use of these products you accept the terms and conditions of all applicable Limited Use Label Licenses. Unless otherwise indicated, these products are for research use only and are not intended for human or animal diagnostic, therapeutic or commercial use.