

## BCIP/NBT Substrate Solution NEL937

### Storage Condition

This product should be stored at 2-8°C.

### Product Description

This phosphatase substrate solution contain 5-bromo-4-chloro-3-indolyl-phosphate (BCIP) at a concentration of 0.21 g/L and nitroblue tetrazolium (NBT) at a concentration of 0.42 g/L in an organic base/Tris buffer. It is provided as 2 x 250 ml of ready-to-use solution, enough to cover ~2000 cm<sup>2</sup> of membrane.

### Performance Characteristics

This ready-to-use substrate solution deposits a permanent dark purple stain on membrane sites bearing phosphatase. Use at room temperature. The amount of substrate solution needed is determined by the size of the membrane being used. Use enough reagent to completely cover the membrane surface or at least 0.25 ml/cm<sup>2</sup> of membrane. It is also useful for in situ hybridization.

### Use of Reagent

1. Incubate at room temperature until a dark purple stain appears at the phosphatase sites.
2. Stop the reaction by rinsing the membrane with reagent quality water to remove excess substrate. Air dry the membrane and store.
3. The point at which the reaction is terminated is determined by the rate of color development. The reaction should be stopped before background color becomes too intense, resulting in insufficient contrast between positive staining and background. To reduce the intensity of the substrate reaction, it is recommended that the phosphatase conjugate and/or antibodies in the assay be further diluted. Dilution of the substrate is not recommended.
4. For long term storage of the stained membrane it is recommended that the dried membrane be overlaid with a clear plastic and stored in the dark.

### Membrane Compatibility

This substrate solution is formulated for use on PolyScreen® PVDF Transfer Membrane, nitrocellulose membranes and the nylon membranes GeneScreen™ Hybridization Transfer Membrane and GeneScreen Plus® Hybridization Transfer Membrane.

### Stability

The product is stable for a minimum of one year at 4°C. A fine black precipitate may develop on storage. The presence of this precipitate does not affect product performance.

### Related Products

NEL300	4CN Plus Chromogenic Substrate	NEF811	Anti-Rabbit IgG (Goat), AP-Labeled
NEL938	DAB Substrate Solution	NEF821	Anti-Mouse IgG (Goat), AP-Labeled
NEL751	Streptavidin-AP Conjugate	NEL761	BLAST® (Chromogenic) For Chromogenic Western Blotting
NEF709	Antifluorescein-AP Conjugate		
NEF801	Anti-Human IgG (Goat), AP-Labeled		