



# **Alkaline Phosphatase Conjugate Substrate Kit**

**Catalog Number  
170-6432**

***BIO-RAD***

**Warning:** This product contains dimethylformamide. See MSDS for important safety information.

## Specifications

---

<b>Contents</b>	<b>AP color reagent A</b> (contains nitroblue tetrazolium in aqueous dimethylformamide [DMF], containing magnesium chloride), 10 ml <b>AP color reagent B</b> (contains 5-bromo-4-chloro-3-indolyl phosphate in DMF), 10 ml <b>AP color development buffer</b> , 25x liquid concentrate, 40 ml
<b>Storage</b>	Store AP color reagents A and B at -20 °C. Store the AP color development buffer solution (both the 25x concentrate and diluted) at 4 °C.
<b>Shelf life</b>	One year for all reagents.

## Directions

1. Color development buffer: Dilute only as much as is needed for the color development reaction, or as much as is practical. The stock solution is a 25x concentrate. Take an appropriate aliquot from the stock solution and dilute to a 1x working solution.

Example dilution scheme:  
To prepare 50 ml of 1x buffer, add 2 ml of 25x concentrate to 48 ml of room temperature deionized water and mix thoroughly. This solution should be used at room temperature. Store any excess 1x solution at 4 °C. Store remaining 25x stock solution at 4 °C.

2. Immediately before use, add 1.0 ml of AP color reagent A and 1.0 ml of AP color reagent B to 100 ml AP color development buffer at RT.
  3. Immerse the membrane in the color development solution. Incubate at RT with gentle agitation until color development is complete.
  4. Stop the development by washing the membrane in dd water for 10 minutes with gentle agitation. Change the water at least once during the 10 minute period to remove residual color development solution.
  5. Air dry the membrane.
- 

This product is for research use only. It is not intended for clinical diagnostic purposes.