

- 2; Immediately before use, add 600 μ l HRP color reagent B to 100 ml of 1x HRP color development buffer at room temperature. Add this solution to 20 ml of HRP color reagent A.
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.



HRP Conjugate Substrate Kit

Catalog Number
170-6431

BIO-RAD

Specifications

Contents	HRP color reagent A (contains 4-chloro-1-naphthol in diethylene glycol), 200 ml HRP color reagent B (contains hydrogen peroxide), 10 ml HRP color development buffer, 100 ml, 10x concentrate
Storage	HRP color reagent A: -20 °C HRP color reagent B: 4 °C HRP color development buffer solution: 4 °C or room temperature. The solution can be frozen. If any salts come out of solution in freezing, warm to room temperature and stir for 10-30 minutes.
Shelf life	One year for HRP color reagents A and B One year for 10x stock HRP color development buffer solution. The 1x working solution should be

made with filtered deionized water. In this condition, the 1x solution can be stored for several months at 4 °C.

- Directions for use**
1. The HRP color development buffer solution is a 10x concentrate. It should be diluted with filtered, deionized water. Dilute to a 1x working solution by adding 1 part HRP color development buffer concentrate to 9 parts filtered, deionized water. For example, to make 100 ml of 1x HRP color development buffer, mix 90 ml of **filtered**, deionized water with 10 ml of 10x HRP color development buffer concentrate. Mix well. Store excess solution at 4 °C.

Since the HRP color development buffer is better stored for longer periods as a concentrate, make as much 1x solution as is practical and needed for the planned experiments.