



DAB ENHANCING SOLUTION

Cat. No. H-2200

Storage: 4°C in the dark

DAB Enhancing Solution is a ready-to-use reagent for toning DAB stained sections. Only visible DAB precipitate is toned, so little amplification is observed with the enhancing reagent.

After developing DAB substrate:

1. Rinse sections in 0.05 M sodium bicarbonate, pH 9.6 for 10 minutes.
2. Carefully blot excess buffer from sections.
3. Cover sections with DAB Enhancing Solution.

NOTE:

Enhancement can take from 5 -20 seconds and should be determined by the investigator. Usually a 10-15 second exposure to the Enhancing Solution is sufficient. The color may vary somewhat depending on the intensity of untoned DAB precipitate. Generally the toned DAB reaction product has a darker reddish brown color. Overexposed tissue will have increased background and may turn pink with time.

4. Quickly rinse sections with water to stop the reaction.
5. Sections can then be counterstained, dehydrated and coverslipped.

Handle with care. Avoid contact with skin.