

DAB Enhancing Solution

Cat. No. H-2200

Storage Refrigerate 2-8 °C

Description DAB Enhancing Solution intensifies the DAB reaction product in stained sections. A ten second exposure to this solution darkens the reaction product.

Components

Product Name	Volume
DAB Enhancing Solution	30 ml

Guidelines for Use

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 and cover sections with DAB Enhancing Solution.
3. Incubate 5-20 seconds. Rinse sections with water.
4. Counterstain sections, if desired, dehydrate, and coverslip.

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.

Handle with care. Avoid contact with skin.