

## 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

<u>Product Name</u>	<u>Volume</u>
DAB Enhancing Solution	30 ml



### Guidelines for Use:

After developing DAB substrate:

Rinse sections in 0.05 M sodium bicarbonate, pH 9.6 for 10 minutes. Carefully blot excess buffer from sections and cover sections with DAB Enhancing Solutions.

**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.

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

Handle with care. Avoid contact with skin.

Detailed product listings, specifications, protocols and additional XXXX.Rev.00