

**VitroView™ DAB Substrate Kit**  
**(for 1000 IHC Tests)**

**Cat. No. VB-6003**

**Introduction**

Diaminobenzidine (DAB) is oxidized by horseradish peroxidase (HRP) to give a dark-brown color. It is a sensitive colorimetric substrate used with HRP in immunohistochemistry (IHC) and immunoblotting applications.

**Kit Components**

1. DAB stock solution (50x): 3 ml
2. Stable H<sub>2</sub>O<sub>2</sub> Solution (50x): 3ml
3. DAB buffer: 120ml

**Storage**

Store at 4°C.

**Application**

Immunohistochemistry: Tested

**Recommended Assay Procedure for Immunohistochemistry**

1. Before DAB staining, add 40 µl of DAB stock solution and 40 µl of Stable H<sub>2</sub>O<sub>2</sub> Solution for every 2 ml of DAB buffer. The mixed DAB substrate solution can be stored for up to 6 hours. Storing the reagent for longer periods of time will cause precipitate formation.
2. Apply the mixed DAB substrate reagent to the tissue section after it has been incubated with the appropriate horseradish peroxidase (HRP) conjugated reagents.
3. Approximately 100-200 µl (2-4 drops) of the mixed DAB substrate reagent should be sufficient to cover the entire tissue section.
4. Incubate the section with the mixed DAB reagent for 2-10 minutes or until the desired color intensity is obtained.
5. Wash the slides in dH<sub>2</sub>O
6. Counterstain

**Warning**

DAB is a possible carcinogen. Please take necessary precautions.

**Precautions**

Handle with care. Avoid contact with eyes, skin and clothing. Do not ingest. Wear gloves.