



## **VitroView™ IHC Hematoxylin Counterstain Kit (300 Tests)**

**Cat. No. VB-6004**

### **Introduction**

This ready-to-used (RTU) product is designed to be used after completion of immunohistochemical staining. The staining color will be light blue in nuclei.

### **Kit Components**

IHC Hematoxylin Counterstain Solution 15ml ×2

### **Storage**

Store at room temperature.

### **Application**

Immunohistochemistry: Tested

### **Recommended Assay Procedure for Immunohistochemistry**

1. After DAB staining, incubate sections with 3 drops of RTU hematoxylin solution for 1-2 minutes
2. Rinse in tap water 2×2 min.
3. Dehydrate through 75% ethanol for 2 min, 95% ethanol for 2 min, and 100% ethanol for 2×3min. Clear in xylene for 2×5min.
4. Coverslip with mounting medium.

### **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.