



## **VitroView™ Antibody Dilution Buffer (100 ml)**

**Cat. No. VB-6002**

### **Introduction**

This antibody dilution buffer is ready-to-use solution for dilution of primary and secondary antibodies in immunoassays including immunofluorescence (IF), Immunohistochemistry (IHC) and Western blot (WB).

### **Kit Components**

BSA (IHC grade, IgG-Free, Protease -free)  
cold fish skin gelatin  
Sodium azide  
0.01M PBS, pH7.2

### **Storage**

Store at 4°C.

### **Application**

IHC, ICC/IF, WB. The usefulness of this product in other applications has not been determined.

### **Note**

1. The sodium azide is an inhibitor of HRP. Do not use it to dilute HRP conjugated antibody.
2. The phosphate in PBS is an inhibitor of Alkaline Phosphatase (AP). Do not use it to dilute AP conjugated antibody.

### **Precautions**

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