

## **EDTA Antigen Retrieval Buffer (1x), 1L**

**Cat. No. VB-6033**

### **Introduction**

The EDTA based solution is designed to break the protein cross-links, therefore unmask the antigens and epitopes in formalin-fixed and paraffin embedded tissue sections, thus enhancing staining intensity of antibodies.

### **Component**

EDTA Antigen Retrieval Buffer (1x), pH 8.0    1L

### **Storage**

Room temperature.

### **Protocol**

1. Deparaffinize sections in 2 changes of xylene, 5 minutes each.
2. Hydrate in 2 changes of 100% ethanol for 2 minutes each, 95% and 75% ethanol for 2 minute each. Then rinse in distilled water.
3. Pre-heat steamer or water bath with staining dish containing 1x EDTA Buffer until temperature reaches 95-100 °C.
4. Immerse slides in the staining dish. Place the lid loosely on the staining dish and incubate for 15 minutes (optimal incubation time should be determined by user).
5. Turn off steamer or water bath and remove the staining dish to room temperature and allow the slides to cool for 20-40 minutes.
6. Rinse sections in PBS Tween 20 for 2x2 min.
7. Proceed to standard immunohistochemistry protocol.

**Note:** Microwave, pressure cooker and autoclave can be used as alternative heating source to replace steamer or water bath.

### **Note**

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