

Citrate Buffer (1x), 1L

Cat. No. VB-6031

Introduction

Citrate buffer is an antigen retrieval buffer for immunostaining.

Component

Citrate Buffer (1x), pH 6.0 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 Citrate 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.