

Antigen Unmasking Solution

Citrate-based (pH 6.0)

Cat. No. H-3300

Storage 2–8 °C

Description Antigen Unmasking Solutions are highly effective at revealing antigens in formalin-fixed, paraffin embedded tissue sections when used in combination with a high temperature treatment procedure.

Components

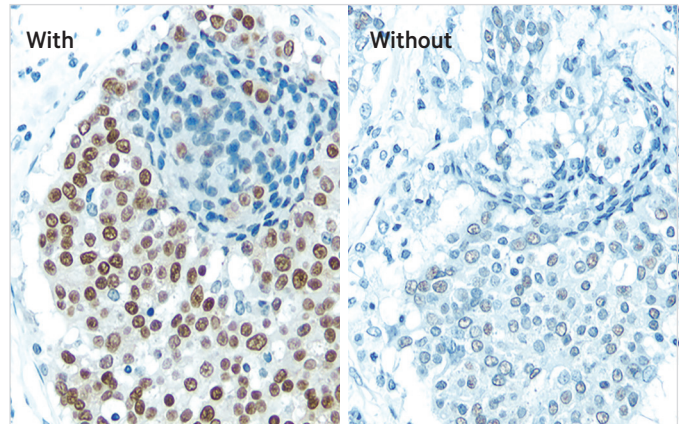
| Product Name | Volume |
|---|--------|
| Antigen Unmasking Solution Citrate-based (pH 6.0) | 250 ml |

Usage Information

Diluted Antigen Unmasking Solutions are designed for use in high temperature unmasking procedures.

The following protocol is for use with a pressure cooker (solution temperature ≈ 120°).

1. Cut and mount sections on slides treated with a tissue section adhesive such as VECTABOND® Reagent*. Position slides into metal staining racks (do not place slides close together; uneven staining may occur).
2. Deparaffinize sections and rehydrate to distilled water.
3. Pour 1600 ml of distilled water into a stainless steel pressure cooker**.
4. Shake well before use, and then add 15 ml of the concentrated stock Antigen Unmasking Solution.
5. Cover, but do not lock lid, and bring solution to a boil.
6. Place slides in pressure cooker ensuring slides are well immersed in diluted Antigen Unmasking Solution. Lock lid.
7. Start timing once the cooker has pressurized, start timing.
8. After one minute, remove the pressure cooker from heat source and run under cold water. DO NOT OPEN LID UNTIL THE INTERNAL PRESSURE HAS BEEN COMPLETELY REDUCED.
9. Open lid, remove slides and place immediately into a tap water bath. DO NOT LET SECTIONS DRY OUT.
10. Wash sections in PBS buffer (pH 7.5) for 5 minutes.
11. Follow standard protocol for immunohistochemical labeling. Many high temperature treatment procedures are compatible with Antigen Unmasking Solutions. Optimize treatment time for individual use.



Breast Carcinoma: with and without Citrate-based (pH 6.0) Antigen Unmasking Solution, Estrogen receptor (er), ImmPRESS® Anti-Rabbit IgG Kit, DAB substrate (brown). Counterstained with Hematoxylin QS (blue).

- * VECTABOND Reagent is a tissue section adhesive that can significantly increase adherence of both frozen and paraffin-embedded tissue sections to glass slides, even under harsh conditions such as those required for high temperature antigen unmasking techniques and in situ hybridization.
- ** Follow the manufacturers instructions and safety recommendations for use.