



Pretreatment Required for using VP-B209 on Paraffin Sections

1. Deparaffinize and rehydrate sections.
2. Incubate with 0.5% hydrogen peroxide in methanol for 10 minutes.
3. Rinse in tap water for 5 minutes.
4. Incubate with trypsin at 37 °C for required time as determined in your laboratory eg 30 minutes.
5. Rinse in tap water for 5 minutes.
6. Preheat in distilled water at 37 °C for 5 minutes.
7. Incubate with 2N HCl at 37 °C for 30 minutes.
8. Incubate with 0.1M di-Sodium tetraborate (Borax) for 10 minutes.
9. Rinse briefly in tap water.
10. Rinse in 50mM Tris-Buffered Saline (pH7.4) for 2 x 5 minutes.
11. Proceed as for standard ABC technique.

Pretreatment Required for using VP-B209 on Frozen Sections

1. Fix in acetone for 10 minutes at 4 °C.
2. Rinse in tap water for 5 minutes.
3. Incubate with 0.5% hydrogen peroxide in methanol for 10 minutes.
4. Rinse in tap water for 5 minutes.
5. Preheat in distilled water at 37 °C for 5 minutes.
6. Incubate with 2N HCl at 37 °C for 30 minutes.
7. Incubate with 0.1M di-Sodium tetraborate (Borax) for 10 minutes.
8. Rinse briefly in tap water.
9. Rinse in 50mM Tris-Buffered Saline (pH7.4) for 2 x 5 minutes.
10. Proceed as for standard ABC technique.