

# Ventana Platform Protocol Summary

## Procedure: U OptiView DAB IHC

### BenchMark ULTRA IHC/ISH

#### IHC Reagents (Ventana Platform):

- Ventana Reaction Buffer (Ventana, Cat# 950-300)
- Optiview Dab IHC Detection Kit (Ventana, Cat# 760-700)
- Ultra CC#1 (Ventana, Cat# 950-224)
- EZ Prep (Ventana, Cat# 950-102)
- LCS (Ventana, Cat# 650-210)
- Antibody Diluent (Dako, Cat# K8006)
- Hematoxylin II (Ventana, Cat# 790-2208)
- Bluing Reagent (Ventana, Cat# 760-2037)

#### Protocol Summary:

1. Paraffin [Selected]
2. Deparaffinization [Selected]
3. Warmup Slide to [72 Deg C] from Medium Temperatures ( Deparaffinization )
4. Cell Conditioning [Selected]
5. Ultra CC1 [Selected]
6. Warmup Slide to [100 Deg C], and Incubate for 4 Minutes ( Cell Conditioner #1 )
7. CC1 8 Min [Selected]
8. CC1 16 Min [Selected]
9. CC1 24 Min [Selected]
10. CC1 32 Min [Selected]
11. CC1 40 Min [Selected]
12. CC1 48 Min [Selected]
13. CC1 56 Min [Selected]
14. CC1 64 Min [Selected]
15. Pre Primary Peroxidase Inhibit. [Selected]
16. Primary Antibody [Selected]
17. Primary Antibody Temperature [Selected]
18. Warmup Slide to [37 Deg C] ( Primary Antibody )
19. Apply Coverslip, One Drop of [PREP KIT 28] ( Antibody ), and Incubate for [0 Hr 16 Min]
20. OptiView HQ Linker [Selected]
21. OptiView HQ Universal Linker [Selected]
22. Apply Coverslip, One Drop of OV HQ UNIV LINKR, and Incubate for [8 Minutes]

23. OptiView HRP Multimer [Selected]
24. HRP Multimer Incubation Time [Selected]
25. Apply Coverslip, One Drop of OV HRP MULTIMER, and Incubate for [8 Minutes]
26. Counterstain [Selected]
27. Apply One Drop of [HEMATOXYLIN II] ( Counterstain ), Apply Coverslip, and Incubate for [8 Minutes]
28. Post Counterstain [Selected]
29. Apply One Drop of [BLUING REAGENT] ( Post Counterstain ), Apply Coverslip, and Incubate for [4 Minutes]

\* one drop is one reagent dispense