

**Labeling Mouse Kidney Sections  
with Qdot® Wheat Germ Agglutinin Conjugate**

## Labeling Mouse Kidney Sections with Qdot® Wheat Germ Agglutinin Conjugate

### PLEASE READ ENTIRE PROTOCOL BEFORE STARTING.

Additional information can be obtained from the product page on our website at [www.invitrogen.com](http://www.invitrogen.com).

### Materials

Acetone fixed mouse kidney sections (Inova Diagnostics, Cat # 508170)  
TBS (Dako, Cat # S3001)  
Tween 20 (Sigma, Cat # P-7949)  
IgG-free BSA (Jackson, Cat # 001-000-162)  
Qdot Wheat Germ Agglutinin Conjugates (Cat # Q12021MP)  
Hoechst 33342 (Molecular Probes, Cat # H3570)  
Polyvinyl Alcohol Mounting Medium with DABCO (Sigma, Cat # 10981)

### Notes:

1. Centrifuge at 5,000-10,000 x g, for 5-10 min. reserving the supernate, prior to using the material.
2. Some PAP pens can quench the signal from the quantum dots. If your protocol requires the use of a PAP pen, we recommend the ImmEdge Hydrophobic Barrier Pen (H-4000) from Vector Labs.

### Procedure

All steps are conducted at room temperature unless specifically indicated. Slides should be placed in a moist chamber for the blocking and incubation steps.

1. Warm the slide to room temperature on bench.
2. Block with 2% BSA/TBS/0.05% Tween-20 for 20 minutes.
3. Incubate with 5 nM Qdot 655 Wheat Germ Agglutinin Conjugate diluted in 2% BSA/TBS at 37 °C for 30 minutes.
4. Wash with TBS for 2 x 3 minutes.
5. Counterstain with 0.5 µg/ml Hoechst/TBS for 5 minutes.
6. Wash with TBS for 2 x 3 minutes.
7. Mount with Polyvinyl Alcohol Mounting Medium/DABCO.