

## NB110-96878 Protocol

### Serum protocol for DNA/RNA Damage Antibody (NB110-96878)

Protocol specific for DNA/RNA Damage Antibody (NB110-96878): [https://www.novusbio.com/products/dna-rna-damage-antibody-15a3\\_nb110-96878](https://www.novusbio.com/products/dna-rna-damage-antibody-15a3_nb110-96878)

#### Tissue Preparation

8-OHdG monoclonal antibody reacts on both 50 um frozen tissue sections and paraffin-embedded sections. Tissue should be dissected fresh and fixed in periodate-lysine-paraformaldehyde (PLP) at 4C over night.

#### PLP

Heat 1 L dH<sub>2</sub>O to 60C.

Add 60 g paraformaldehyde.

Add 33 g dibasic NaPO<sub>4</sub>.

Cool to room temperature in a cold water bath.

Add 9 g monobasic NaPO<sub>4</sub>.

Add 6.45 g Na-m-periodate.

Add 41.1 g lysine (HCl salt).

Filter and dilute to 3 L with dH<sub>2</sub>O.

Adjust pH to 7.6 with 1.0 N NaOH approx. (20-30 ml).

Tissue prepared for frozen sectioning must be cryoprotected in a 20% glycerol-2% DMSO solution in phosphate buffer for 24-48 hours. Tissue will sink to the bottom of container when fully penetrated. This will eliminate freezing artifact from cutting.

Glycerol-DMSO (for 3 L)

2.4 L 0.1 M phosphate buffer

600 ml glycerol

60 ml DMSO

0.1 M Phosphate Buffer, pH 7.4 (for 1 L)

1 L dH<sub>2</sub>O

11 g dibasic NaPO<sub>4</sub>

3 g monobasic NaPO<sub>4</sub>

After frozen sectioning, tissue should be stored in phosphate buffer with 0.08% sodium azide.

#### Staining Sections By DAB Procedure

Paraffin-embedded sections must be deparaffinized by sequential immersion in the following for 3 minutes each: xylene (twice), absolute ethanol (twice). Agitate gently in each solution. Proceed with the following procedure.

1. Pretreat sections with a methanol-peroxide solution to eliminate endogenous peroxidases.

Methanol-Peroxide

100 ml absolute methanol

1 ml 33% H<sub>2</sub>O<sub>2</sub>

Incubate sections in methanol-peroxide solution for 30 minutes, room temperature.

2. Wash sections 3 times for 10 minutes each in 0.1 M phosphate buffered saline (PBS)

PBS, pH 7.4 (for 1 L)  
1 L dH<sub>2</sub>O  
11 g dibasic NaPO<sub>4</sub>  
3 g monobasic NaPO<sub>4</sub>  
8.5 g NaCl

3. Incubate sections for 1 hour in 10% normal goat serum in PBS.

4. Incubate sections in the primary antibody for 18-24 hours at room temperature. Depending on the nature of the sample, a shorter incubation time may be used. It is recommended that a concentration range of 1-10 ug/ml be evaluated in order to determine the optimal concentration for each type of tissue sample. Dilute antibody in PBS containing 0.3% Triton X-100, 0.08% sodium azide and 2% normal goat serum.

NOTE: A humidified chamber is necessary when staining paraffin sections. Slides should be placed flat and primary antibody applied over the section, covering it completely.

5. Rinse sections 3 times for 10 minutes each in PBS.

6. Incubate for 3 hours with peroxidase-conjugated goat anti-mouse IgG (Boehringer-Mannheim, Indianapolis, IN) diluted 1:300 in PBS with 2% normal goat serum.

7. Rinse sections 3 times for 10 minutes each in PBS.

8. Incubate sections for 5-10 minutes in a solution of 0.5 mg/ml 3,3' diamino-benzidine tetrahydrochloride (DAB, Sigma, St. Louis, MO) and 0.005% hydrogen peroxide in 0.05 M tris HCl buffer, pH 7.6 plus imidazole (10 ml/110 ml Tris buffer).

50 mM Tris Buffer, pH 7.6  
1 L dH<sub>2</sub>O  
6 g Trizma base  
3 ml concentrated HCl (37%)

Sodium Imidazole  
100 ml 0.1 M phosphate buffer  
0.7 g sodium imidazole

9. Rinse sections 3 times for 10 minutes each in PBS.

10. Mount free-floating sections on subbed slides and air dry.

Subbing Solution  
500 ml dH<sub>2</sub>O

2.5 g gelatin

0.25 g chromium potassium sulfate

Heat to 60°C. Filter and proceed to coat slides. Once slides are air dried, sections can be mounted.

11. Dehydrate mounted/paraffin sections by sequential immersion in the following for 3 minutes each: 70% ethanol, 95% ethanol, absolute ethanol, xylene. Agitate gently in each solution.

12. Apply coverslip with Permount in a chemical fume hood