

NB21-1002 Protocol

Protocol specific for H3R2me2a antibody (NB21-1002)

Histone H3 [Asym-dimethyl Arg2] Antibody: https://www.novusbio.com/products/histone-h3-antibody_nb21-1002
Immunocytochemistry Protocol

Culture cells to appropriate density in 35 mm culture dishes or 6-well plates.

1. Remove culture medium and add 10% formalin to the dish. Fix at room temperature for 30 minutes.
 2. Remove the formalin and add ice cold methanol. Incubate for 5-10 minutes.
 3. Remove methanol and add washing solution (i.e. PBS). Be sure to not let the specimen dry out. Wash three times for 10 minutes.
 4. To block nonspecific antibody binding incubate in 10% normal goat serum from 1 hour to overnight at room temperature.
 5. Add primary antibody at appropriate dilution and incubate at room temperature from 2 hours to overnight at room temperature.
 6. Remove primary antibody and replace with washing solution. Wash three times for 10 minutes.
 7. Add secondary antibody at appropriate dilution. Incubate for 1 hour at room temperature.
 8. Remove antibody and replace with wash solution, then wash for 10 minutes. Add Hoechst 33258 to wash solution at 1:25,000 and incubate for 10 minutes. Wash a third time for 10 minutes.
 9. Cells can be viewed directly after washing. The plates can also be stored in PBS containing Azide covered in Parafilm (TM). Cells can also be cover-slipped using Fluoromount, with appropriate sealing.
- *The above information is only intended as a guide. The researcher should determine what protocol best meets their needs. Please follow safe laboratory procedures.