



Orders: orders@novusbio.com

Support: technical@novusbio.com

Web: www.novusbio.com

Protocols, Publications, Related Products, Reviews and more:

www.novusbio.com/NBP2-29865

NBP2-29865 Protocol

Serum protocol for MATH2/NEUROD6 Antibody (NBP2-29865)

Protocol specific for MATH2/NEUROD6 Antibody (NBP2-29865): https://www.novusbio.com/products/math2-neurod6-antibody_nbp2-29865

To reconstitute the antibody, centrifuge the antibody vial at moderate speed (5,000 rpm) for 5 minutes to pellet the precipitated antibody product. Carefully remove the ammonium sulfate/PBS buffer solution and discard. It is not necessary to remove all of the ammonium sulfate/PBS solution: 10 uL of residual ammonium sulfate solution will not effect the resuspension of the antibody. Do not let the protein pellet dry, as severe loss of antibody reactivity can occur.

Resuspend the antibody pellet in any suitable biological buffer, standard PBS or TBS (pH 7.3-7.5) are typical. Volumes required are not critical but it is suggested that the final antibody concentration be between 0.1 mg/mL and 1.0 mg/mL. For example, to achieve a 1 mg/mL concentration with 50 ug of precipitated antibody, the amount of buffer needed would be 50 uL.

Carefully add the liquid buffer to the pellet. DO NOT VORTEX. Mix by gentle stirring with a wide pipet tip or gentle finger-tapping. Let the precipitated antibody rehydrate for 1 hour at 4-25C degrees prior to use. Small particles of precipitated antibody that fail to resuspend are normal. Vials are overfilled to compensate for any losses.