

Reconstitution Instructions for NBP3-18318

Dried beta Amyloid 42 peptide has been pre-treated with hexafluoroisopropanol (HFIP) to disrupt any fibrils formed after peptide synthesis. Novus Biologicals recommends re-suspended peptide is kept on ice and used immediately to avoid aggregation. While several methods of monomer resuspension are detailed in the literature, Novus Biologicals recommends the following protocol for re-suspending 100 µg of peptide into a 1 mg/mL solution:

1. Bring dried peptide to room temperature for 10 minutes.
2. Add 7.5 µL cold filter sterilized 1% Ammonium hydroxide (NH₄OH) to the peptide aliquot, mix well by pipetting up and down. Take care to scrape inside lower walls of tube to ensure complete re-suspension of the film.
3. Add 92.5 µL of cold filter sterilized PBS. *Optional - Centrifuge monomer 14,000 x g at 4°C for approximately 5-10 minutes and keep the supernatant to remove any material that was not fully re-suspended.*
4. Keep monomer on ice and use immediately.

While it is not recommended, depending on your application small aliquots (≤ 50 µL) can be frozen and kept for up to 2 weeks at -80°C. However, aggregate should be removed by centrifugation at 4°C upon thawing.