

Is there a stop solution or fixative that can be used to stop a reaction prior to analysis on the Luminex analyzer?

We have used PBS-0.2% Formaldehyde (150 mM Sodium Chloride, 10 mM Sodium Phosphate, 0.2% v/v Formaldehyde, pH 7.4) as a stop solution to fix reactions before reading on the Luminex analyzer. **Note:** It may be necessary to titrate the amount of formaldehyde for the specific application.

A filter plate is required for this procedure. Assays not performed in a filter plate must be transferred prior to fixation.

1. After incubation with the reporter fluorophore is completed, transfer the reactions to a pre-wetted filter plate if necessary:
 - a. Pre-wet the appropriate wells of a 1.2 μm Millipore filter plate with 100 μL /well of PBS-1% BSA and aspirate by vacuum manifold.
 - b. Add 50 μL of PBS-1% BSA to the appropriate wells of the filter plate.
 - c. Transfer the contents of the assay plate to the filter plate.
2. Aspirate the supernatant by vacuum manifold.
3. Wash each well twice with 100 μL PBS-1% BSA and aspirate by vacuum manifold.
4. Add 50 μL of PBS, pH 7.4 to each well.
5. Add 50 μL of PBS-0.2% Formaldehyde to each well.
6. Resuspend the microspheres by gently pipetting up and down several times using a multi-channel pipettor.