

Preparation of Antigen Standard

Preparation instructions for Kits containing more than 5 different antigen standard sets

Different antigen standard set vials can be reconstituted simultaneously as long the volume of sample type-specific standard buffer is at least 50 μL per vial and equals 250 μL in total. For your convenience an example schema is included below.

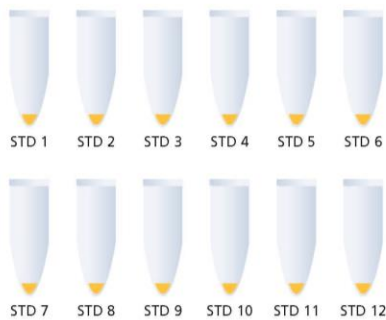
Step 1. Reconstitution and pooling of Standards

- Centrifuge all different antigen standard set vials at 2000 $\times g$ for 10 sec.
- Add 50 μL sample type-specific standard buffer into 5 different standard set vials. For serum or plasma samples, use Universal Assay Buffer and for cell culture supernatant samples use the cell culture media that was used to culture the cells.
- Gently vortex the vial(s) for 30 seconds and centrifuge at 2000 $\times g$ for 10 seconds to collect contents at the bottom of the vial(s).
- Incubate on ice for 10 min to ensure complete reconstitution.
- Transfer the entire contents of each vial into different standard set vial and repeat steps C-E until all different antigen standard vials are reconstituted.
- Pool entire contents of each vial into one of the vials.
- Gently vortex the vial for 30 seconds and centrifuge at 2000 $\times g$ for 10 seconds to collect contents at the bottom of the vial.

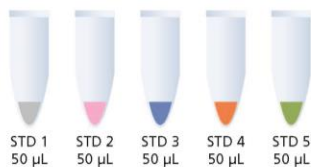
After reconstitution is finished proceed to "Step 2. Prepare 4-Fold Serial Dilution" under Antigen Standard in ProcartaPlex user manual.

Example for reconstitution of 12 different antigen standard set vial

1. Centrifuge all different antigen standard vials



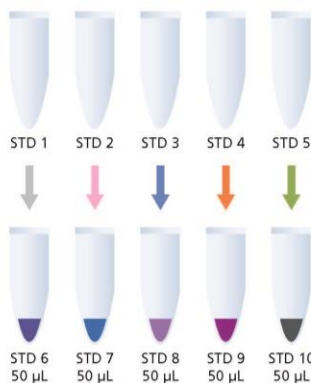
2. Add sample type-specific standard buffer



3. Vortex and centrifuge

4. Incubate on ice

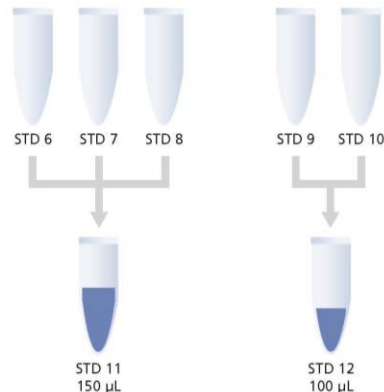
5. Transfer the whole content into different standard vials



6. Vortex and centrifuge

7. Incubate on ice

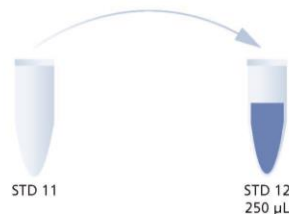
8. Transfer the whole content into different standard vials



9. Vortex and centrifuge

10. Incubate on ice

11. Transfer the whole content into one of the vials



12. Vortex