

Preparation of Antigen Standard

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

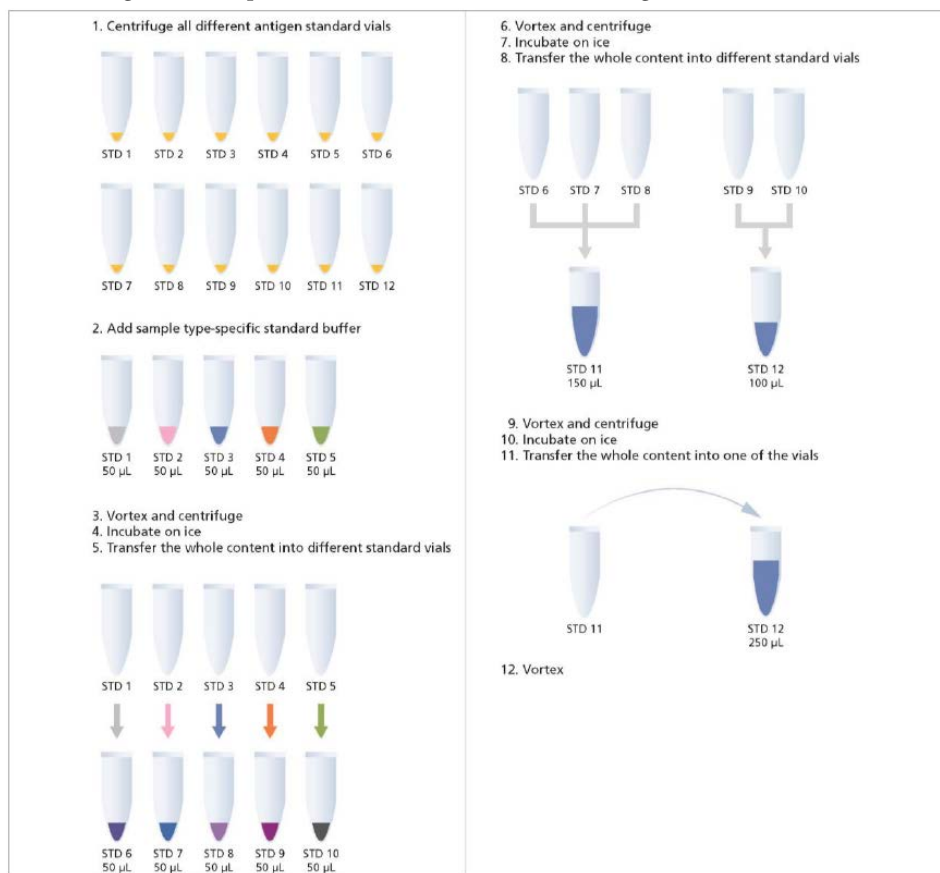
Different antigen standard set vials can be reconstituted simultaneously as long as the volume of sample type-specific standard buffer is at least 50 μL per vial and equals 250 μL in total. For your convenience, an example schema is included below.

Step 1. Reconstitution and pooling of Standards

1. Centrifuge all different antigen standard set vials at $2,000 \times g$ for 10 seconds.
2. Add 50 μL of 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 used to culture the cells.
3. Gently vortex the vial(s) for 30 seconds and centrifuge at $2,000 \times g$ for 10 seconds to collect contents at the bottom of the vial(s).
4. Incubate on ice for 10 minutes to ensure complete reconstitution.
5. Transfer the entire contents of each vial into different standard set vial and repeat steps 3-5 until all different antigen standard vials are reconstituted.
6. Pool entire contents of each vial into one of the vials. Final total volume should equal 250 μL .
7. Gently vortex the vial for 30 seconds and centrifuge at $2,000 \times g$ for 10 seconds to collect contents at the bottom of the vial.

After reconstitution is completed, proceed to "Step 2. Prepare 4-fold serial dilution" under Antigen Standard in the ProcartaPlex™ user manual.








Figure 1. Example for reconstitution of 12 different antigen standard set vials.



Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at www.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have any questions, please contact Life Technologies at www.thermofisher.com/support.

Product label explanation of symbols and warnings

 REF	Catalog Number	 LOT	Batch code		Temperature limitation		Use by		Manufacturer		Consult instructions for use		Caution, consult accompanying documents
--------------------------------------------------------------------------------------	----------------	---------------------------------------------------------------------------------------	------------	-----------------------------------------------------------------------------------	------------------------	-----------------------------------------------------------------------------------	--------	-----------------------------------------------------------------------------------	--------------	-------------------------------------------------------------------------------------	------------------------------	-------------------------------------------------------------------------------------	-----------------------------------------

Manufacturer's address: Bender MedSystems GmbH | Campus Vienna Biocenter 2 | 1030 Vienna, Austria

The information in this guide is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Important Licensing Information: These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2018 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.