

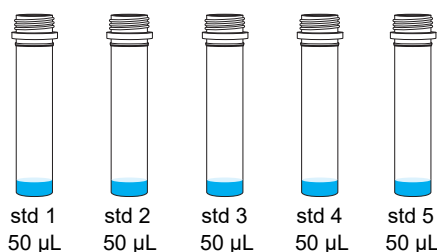
Preparation of a working standard for ProcartaPlex™ Mix & Match panels with more than 5 standards

Pub. No. MAN0017832 Rev. B.0

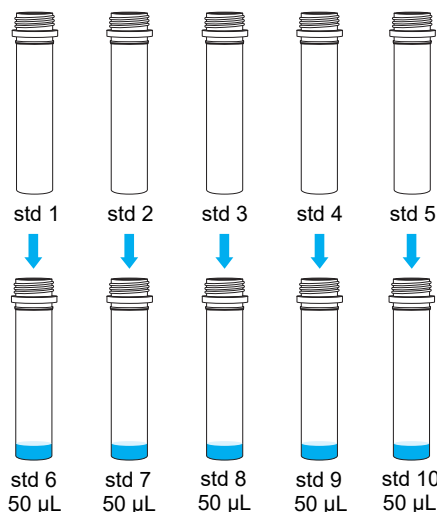
WARNING! Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves. Safety Data Sheets (SDSs) are available from thermofisher.com/support.

This protocol demonstrates the procedure for reconstituting and pooling 12 antigen standard vials, but can be modified for any number of standards greater than 5. Each vial needs to be reconstituted in at least 50 µL and the total volume at the end will be 250 µL. A video demonstrating the procedure by mixing 6 antigen standard vials is available at thermofisher.com/multivial-antigen-prep

1. Remove one of each standard stock vial. Centrifuge each vial at 2,000 x g for 10 seconds.
2. Choose the first 5 standard stock vials (std 1-5 below) and open carefully on the lab bench. Depending on the sample type, add 50 µL of either 1X UAB or cell culture medium. Vortex all 5 vials at high speed for 30 seconds.

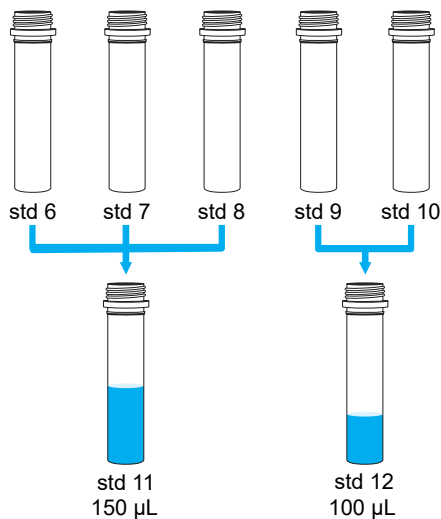


3. Centrifuge at 2,000 x g for 10 seconds to collect contents at the bottom of the vial.
4. Incubate on ice for 10 minutes to ensure complete reconstitution.
5. Transfer 50 µL from each reconstituted vial into the next 5 standard stock vials (std 6-10 below) and vortex the vials at high speed for 30 seconds.

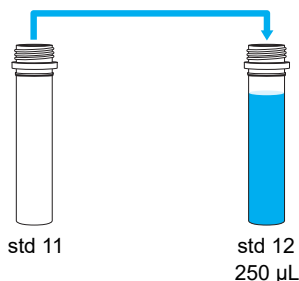


6. Centrifuge at 2,000 x g for 10 seconds to collect contents at the bottom of the vial.
7. Incubate on ice for 10 minutes to ensure complete reconstitution.

8. Transfer 50 μL of each of the 5 reconstituted standard vials into the remaining 2 standard stock vials (std 11-12 below) and vortex the vials at high speed for 30 seconds.



9. Centrifuge at $2,000 \times g$ for 10 seconds to collect contents at the bottom of the vials.
10. Incubate on ice for 10 minutes to ensure complete reconstitution.
11. Pool the contents of the 2 vials (std 11-12 below) into a single vial so the final volume should be 250 μL .



12. Vortex the working antigen standard vial at high speed and then centrifuge at $2,000 \times g$ for 10 seconds to collect contents at the bottom of the vial.



Bender MedSystems GmbH | Campus Vienna Biocenter 2 | 1030 Vienna, Austria
 For descriptions of symbols on product labels or product documents, go to [thermofisher.com/symbols-definition](https://www.thermofisher.com/symbols-definition).

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

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. 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.

Revision history: Pub. No. MAN0017832

Revision	Date	Description
B.0	7 July 2021	Updated the procedure to align with user manuals for ProcartaPlex™ Mix & Match panels.
A.0	11 June 2018	New manual.

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.

©2021 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.