

August 2012

Commercial Process Controls and Re-Gating Samples after a Run with BD Multiset™ Software

How do we analyze and gate commercial process controls with BD Multiset software?

Commercial process controls frequently have dimmer than normal staining. For this reason, College of American Pathologists (CAP) specimens might need to be adjusted more frequently, since these cells are fixed, as are the process controls. In the product insert supplied with the CAP samples, visual inspection and manual gating are recommended when running the proficiency-testing specimens.

It is permissible to adjust the SSC and threshold on the samples to give the best separation between lymphs and monos. It is important to optimize the acquisition to give the software the best opportunity to analyze the data correctly.

Since the attractors are defined based on staining patterns observed with fresh whole human blood, the CD3⁺ attractor might need to be adjusted on all tubes of a multitube panel. Verify that each subset is properly classified. See Figures 1a, 1b, and 1c. Note the inadequate resolution between the CD3⁺ and CD3⁻ lymphocytes in Figure 1b.

The first step that the software takes is to identify the bead population using the CD3/8/45/4 (if BD Trucount™ tubes are used). It is important not to include any cells from the CD3⁺ and CD8⁺ double positives, which can encroach on this population. It is also important to make sure that all of the beads are included within the bead gate. This is a critical gate because if the number of beads is falsely elevated it will make the calculation for absolute counts incorrect. Each entire distinct population should be represented by one unique color.

The second step that the software takes is to place the CD3 attractors, both CD3⁺ and CD3⁻. Once an event is classified by a primary attractor (bead or CD3⁺ or CD3⁻), it cannot be classified by another, even if another attractor overlaps it.

Review the Expert Gate. If events are included that should be eliminated, or if events were missed that should be included, then regate. Sometimes regating is necessary to get the cleanest lymphocyte gate. The operator can always return to the automatic expert gate.

Review the CD3⁺ gate (CD3 vs SSC). Remember that the CD3⁺ population can be dimmer with fixed samples, but do not adjust the gate to encroach into the CD3⁻ population as the gate is shown doing in Fig. 1b.



August 2012

Commercial Process Controls and Re-Gating Samples after a Run with BD Multiset™ Software

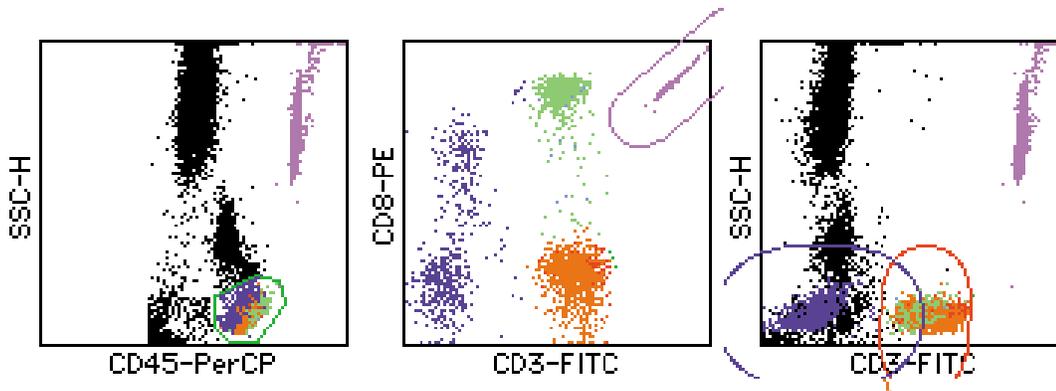


Figure 1a. Fresh whole blood sample showing adequate resolution between the CD3⁺ and CD3⁻ lymphocytes

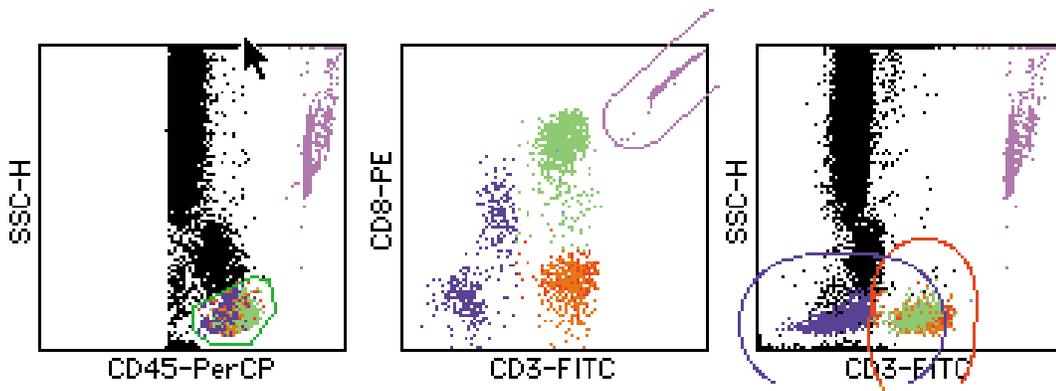


Figure 1b. Preserved process control *before* the CD3⁺ attractor correction

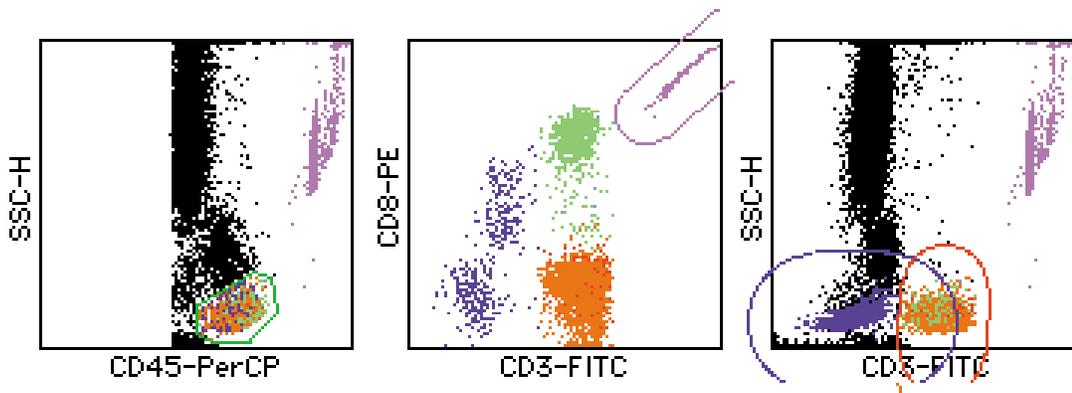


Figure 1c. Preserved process control *after* the CD3⁺ attractor correction

August 2012

Commercial Process Controls and Re-Gating Samples after a Run with BD Multiset™ Software

Finally, look at the subset attractors in the CD8 PE vs CD4 APC plot. Verify that the CD4⁺ and CD8⁺ attractors are in the correct position: the CD3⁺,CD4⁺ attractor followed by the CD3⁺,CD8⁺ attractor and finally the CD3⁺,CD4⁺,CD8⁺ attractor. It is OK for them to overlap, as long as the CD3⁺,CD4⁺ attractor covers all the CD4⁺ cells, and the CD3⁺,CD8⁺ attractor covers all the CD8⁺ cells. The CD3⁺,CD4⁺,CD8⁺ attractor will likely overlap both of the other attractors in this plot, but the true CD4 and CD8 cells have already been captured by their appropriate attractors before this attractor was placed. The CD3⁺,CD4⁺,CD8⁺ attractor is supposed to capture all events that are dual-positive for CD4 and CD8. If any of the attractors needs to be adjusted to enclose their appropriate populations of interest, then do so.

For the 3/16 + 56/45/19 tube, again the bead attractor is reviewed and adjusted if necessary. Check the expert gate and adjust if necessary. Check and adjust the CD3⁺ attractor *and* the CD3⁻ attractor. In this tube, the CD3⁻ cells are the ones that will be considered to be B or NK cells. Next, review and adjust the CD19⁺ attractor (B cells) and then the CD16⁺,CD56⁺ attractor (NK cells).

Observe any QC messages. The lymphosum should fall within a range of 95 to 105. If not, this is a clue to troubleshoot the results. Again, adjustment of the Expert Gate and attractor positions will most likely be required on fixed samples such as process controls and CAP survey samples.

