

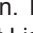
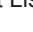


## Time to Analyze!

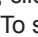

The Analyze tab is used to display multiple plots and samples in any combination, supporting easy analysis.

### Make a New Plot

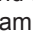

1. Click on an open plot grid, and then click on the Histogram , Dot Plot , Density Plot , or Overlay Histogram  button. The new plot displays in the Plot List.
2. Select a plot and click on a sample in the 96-well grid to view data.

*Tip! The Analyze tab allows simultaneous display of multiple samples.*

### Hide or Delete a Plot

1. To hide a plot from view, click on the Minimize  button. To show it again, click on the plot name in the plot list.
2. To delete a plot from the plot list, click on the Delete  button.

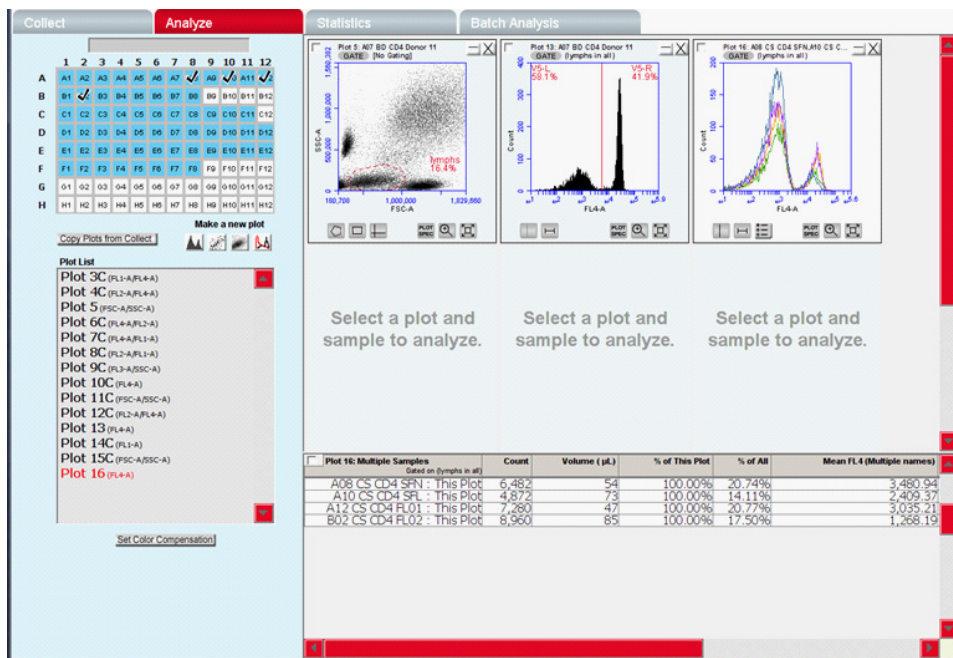
### Make an Overlay Histogram

1. Click on an open plot grid and then click on the Overlay Histogram  button.
2. Select up to 96 samples in the well grid.
3. Click on the Legend  button to view or change the sample color assignments.

*Tip! You can drag plots from C6 Software to other applications.*

### Print Multiple Plots

1. Click on the checkbox in the upper left hand corner of the plot or the statistics table to be printed.
2. Select the **Print Selected Items** option in the File menu.



### Copy a Plot from the Collect Tab

1. Click on the **Copy Plots from Collect** button.
2. In the **Copy Plots** dialog, select the plots that you want to display in the Analyze tab. The plot names display in the **Plot List** followed by **-C** to indicate copies.
3. Click on an open plot grid, select the plot that you want to display, and then click on a sample in the well grid. Gating information from Collect does not copy to Analyze.

### Set VirtualGain™

**VirtualGain** is used to standardize the position of a given peak from sample to sample.

1. On a histogram plot, click on the x-axis parameter name and then click on the **VirtualGain** option from the drop-down menu.
2. In the **VirtualGain** dialog, select the sample that you want to align, move the markers to define the peak, and then click on the **Preview** button. When **VirtualGain** is applied as desired, click on the **Apply** button. A black asterisk \* indicates Gained data.