

## Developing Your Statistics

The Statistics tab is used to display statistics for some or all of your collected samples. You will need at least one sample and a plot to view statistics.

- A** **Preview** shows a preview image of the selected plot/sample combination.
- B** **Display Plot Preview** lists all plots created on the Collect and Analyze tabs.
- C** **Statistics Column Selector** displays all available plot names, their gates, and associated statistics columns.
- D** **Sample Selector** displays all samples collected or imported.
- E** **Master Statistics Table** shows a custom table of statistics selected from Column and Sample Selectors.

### Build a Statistics Table

1. Include a statistic in your table by clicking the corresponding check box in the **Statistics Column Selector** area.
2. Uncheck a column to remove it from the table.

### Preview a Plot or Sample

1. Use Display Plot Preview to select a plot to be previewed.
2. Use **Preview** column of the **Sample Selector** to select a sample to be previewed.

*Tip! Previews are for viewing purposes only and cannot be edited from the Statistics tab.*

### Select Samples

1. Select the sample to be included by clicking on the check boxes from the 'Add to Table' column.
2. The sample names are displayed and the statistics are added to the **Master Statistics Table**.
3. Uncheck samples to remove them from the **Master Statistics Table**.

**Sample Selector**  
Add rows to your master statistics table by selecting samples.

Preview	Add to Table	Sample Name
<input type="checkbox"/>	<input checked="" type="checkbox"/>	HPB
<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	HPB Auto, CD11b,C...
<input type="checkbox"/>	<input checked="" type="checkbox"/>	HPB CD3-F CD4, CD...

### Master Statistics Table

1. The Master Statistics Table is saved as part of the workspace by saving the BD Accuri C6 Software file.
2. You can also copy and paste statistics into spreadsheet applications for calculations or printing purposes.

*Tip! This is a convenient way to record validation bead data into a spreadsheet for tracking instrument performance.*