

AMPLIFICATION CFX Manager™ Software Plate Quick Guide

Run Setup Plate Tab

The Plate tab displays a preview of the plate loaded in the Run Setup window (Figure 1).

Click **Create New** to open the Plate Editor to create a new plate.

Click **Select Existing** to launch the file browser to load a plate file to use in a run or to edit.

Use the Express Load dropdown menu to directly load a plate file to use in a run or to edit.

Click **Edit Selected** to open the Plate Editor to edit the well contents of the selected plate.

Click the **Start Run** tab to proceed and start a run with the currently loaded plate.

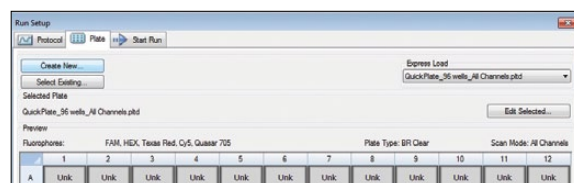


Fig. 1. Plate tab in the Run Setup window. Load an existing plate or create a new plate for a run.

Plate Editor

The Plate Editor is used to create a new plate or edit an existing one (Figure 2).

1. Use the Scan Mode dropdown menu in the Plate Editor toolbar to designate the data acquisition mode to be used during the run.
2. Click **Select Fluorophores** to indicate the fluorophores that will be used in the run.
3. Within the plate diagram, select the wells to load.
4. Choose the Sample Type from the dropdown menu.
5. Click the appropriate checkbox(es) to load the fluorophore(s) in the selected wells.
6. Type the Target Name for each fluorophore (required for gene expression analysis) and press **Enter**, or choose one from the dropdown menu.
7. Type the Sample Name (required for gene expression analysis) and press **Enter**, or choose one from the dropdown menu.
8. To enter the Biological Set Name, check **Biological Set** in the View box at the bottom of the data analysis window.
9. For gene expression analysis, click **Experiment Settings** to assign reference targets and a control sample.

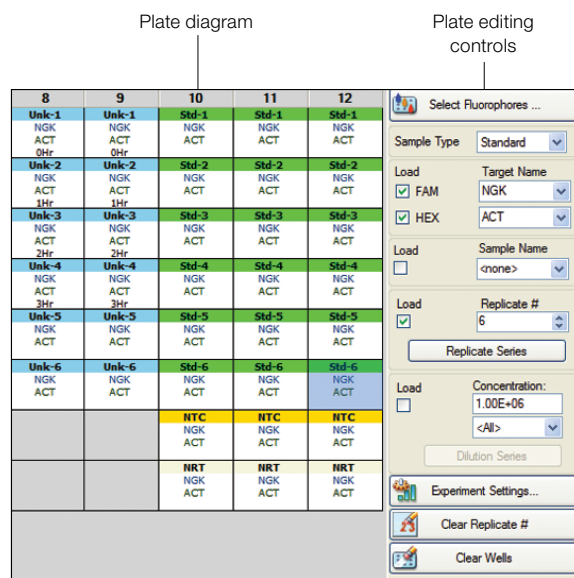


Fig. 2. Plate Editor. In the plate diagram, select the wells you want to load. Use the plate editing controls to enter or edit well contents.

Entering Replicate Numbers

To designate a set of replicate wells, highlight the wells and type or choose a replicate number in the Replicate # box in the plate editing controls (Figure 3). Alternatively, to assign replicate numbers to several well subsets at once:

1. Select wells in the plate diagram and click **Replicate Series**. The Replicate Series editing window opens (Figure 3).
2. Enter the Replicate Group Size and the Starting Replicate #.
3. Indicate whether replicates are loaded horizontally or vertically.
4. Click **Apply** to enter the replicate numbers.

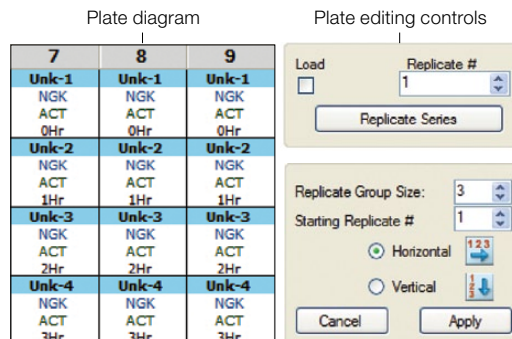


Fig. 3. Replicate Series editing window opens in the plate editing controls.

Creating a Standard Curve

To enter the starting target concentration of one standard, select the wells loaded with Sample Type/Standard, enter a value under Concentration in the plate editing controls (Figure 4), designate All or a specific fluorophore, and then click the **Load** checkbox. Alternatively, to enter concentrations for the entire standard curve series at once:

1. Select the wells that have also been assigned consecutive replicate numbers and click **Dilution Series**. The Dilution Series window opens (Figure 4).
2. Enter the Starting Concentration of the dilution series.
3. Enter the numbers of the first and last replicates in the series.
4. Enter the Dilution Factor and indicate whether the dilution is increasing or decreasing (that is, whether the value entered in Step 2 is the lowest or highest concentration).
5. Click **Apply** to assign the dilution series.

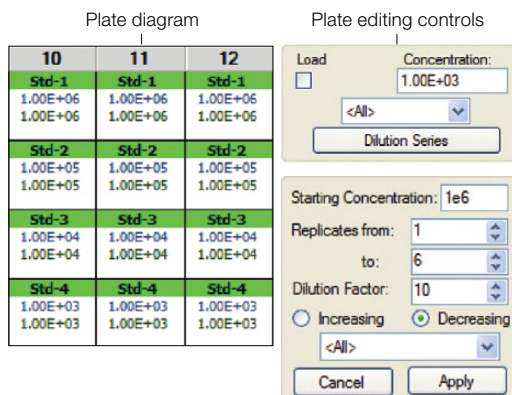


Fig. 4. Dilution Series window for creating a standard curve opens in the plate editing controls.

Creating Well Groups

To create well groups that are analyzed independently:

1. Click the **Well Groups** button in the Plate Editor toolbar. The Well Groups Manager window opens (Figure 5).
2. Click **Add** to create a new group.
3. In the plate diagram, select the wells that will constitute the well group.
4. Click **OK** to return to the Plate Editor window.

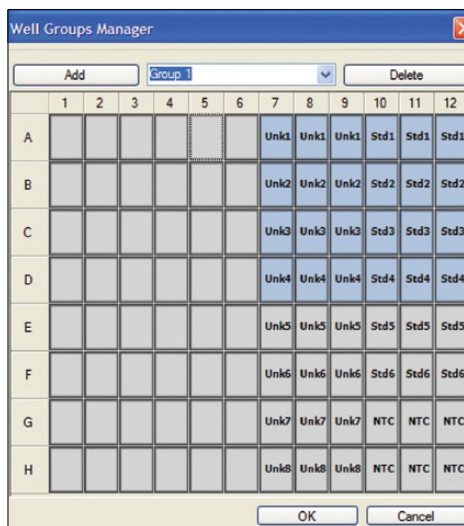


Fig. 5. Well Groups Manager window.

Bio-Rad's real-time thermal cyclers are covered by one or more of the following U.S. patents or their foreign counterparts owned by Eppendorf AG: U.S. Patent Numbers 6,767,512 and 7,074,367.

For more information, visit www.bio-rad.com/web/ampSW96plate.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com USA 800 424 6723 Australia 61 2 9914 2800 Austria 01 877 89 01 Belgium 09 385 55 11 Brazil 55 11 5044 5699
Canada 905 364 3435 China 86 21 6169 8500 Czech Republic 420 241 430 532 Denmark 44 52 10 00 Finland 09 804 22 00
France 01 47 95 69 65 Germany 089 31 884 0 Greece 30 210 9532 220 Hong Kong 852 2789 3300 Hungary 36 1 459 6100 India 91 124 4029300
Israel 03 963 6050 Italy 39 02 216091 Japan 03 6361 7000 Korea 82 2 3473 4460 Mexico 52 555 488 7670 The Netherlands 0318 540666
New Zealand 64 9 415 2280 Norway 23 38 41 30 Poland 48 22 331 99 99 Portugal 351 21 472 7700 Russia 7 495 721 14 04
Singapore 65 6415 3188 South Africa 27 861 246 723 Spain 34 91 590 5200 Sweden 08 555 12700 Switzerland 026 674 55 05
Taiwan 886 2 2578 7189 Thailand 800 88 22 88 United Kingdom 020 8328 2000