

AMPLIFICATION

CFX Manager™ Software Gene Expression Analysis Quick Guide

With stringently qualified controls, you can use the Gene Expression Module of CFX Manager software to evaluate relative differences in a target's concentration between samples. The most common use of this application is for evaluating the target concentration in cDNA samples to infer steady-state mRNA levels. Typically, the message level(s) of one or more reference genes is used to normalize the values of a gene of interest. Reference genes correct for loading differences or other sampling variations present in each sample. The expression level of reference genes should not vary across the biological conditions being studied.

Gene Expression Analysis Setup

The example wells shown in Figure 1 contain a single fluorophore loaded in the wells with four replicate groups containing two different target names and two different sample names.

Assigning Reference Genes

1. Click **Experiment Settings** in the Plate Editor or the Gene Expression Module. The Experiment Settings window opens (Figure 2).
2. Select the **Targets** tab.
3. Click the appropriate checkbox(es) to indicate the target(s) that will be used as references.
4. Click on the **Show Analysis Settings** checkbox in order to use a reaction efficiency other than 100% for any target.
5. The Auto Efficiency checkbox is checked by default. If a standard curve has been included in the run, then the efficiency calculated from the standard curve will be used in the calculations. Alternatively, type a reaction efficiency value into the cell corresponding to the appropriate target to use this efficiency in the calculations.

Assigning Control Samples

1. In the Experiment Settings window, select the **Samples** tab.

	1	2	3	4
A	Unk-1 Target 1 Sample 1	Unk-2 Target 1 Sample 2	Unk-3 Target 2 Sample 1	Unk-4 Target 2 Sample 2
B	Unk-1 Target 1 Sample 1	Unk-2 Target 1 Sample 2	Unk-3 Target 2 Sample 1	Unk-4 Target 2 Sample 2
C	Unk-1 Target 1 Sample 1	Unk-2 Target 1 Sample 2	Unk-3 Target 2 Sample 1	Unk-4 Target 2 Sample 2

Fig. 1. Representative wells for gene expression.

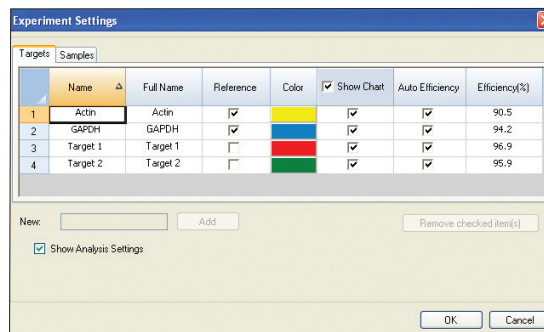


Fig. 2. Assigning reference target(s).

- Click the appropriate checkbox to indicate the sample that will be used as the control in the calculations (Figure 3). The control sample is assigned a value of 1 for every gene, while all other samples are presented with values relative to the control.

Gene Expression Analysis

Selecting the Analysis Mode

- Select the analysis mode from the Mode drop-down menu (Figure 4):
 - Normalized expression $\Delta\Delta C_q$ — relative quantity of gene(s) of interest is normalized to relative quantity of the reference gene(s) across samples.
 - Relative quantity ΔC_q — relative quantity of gene(s) of interest is not normalized; the results express the gene concentration relative to other samples in the experiment.

Graphing Options

- Select from the Graph Data drop-down menu to display graph data relative to control (originating at 1) or relative to zero. In log scale, the Graph Data option defaults to relative to control.
- Select from the X-axis drop-down menu to display Sample or Target on the x-axis.
- Select from the Y-axis drop-down menu to use Linear, Log 2, or Log 10 as the y-axis scale.
- When needed, select from the Scaling drop-down menu — Highest, Lowest, or Unscaled.
- Select from the Error Type drop-down menu to display Standard Error of the Mean or Standard Deviation.
- Choose the Sort option on the right-click graph menu to customize the presentation order on the x-axis.
- Right click on the graph to copy and paste it into other applications or to save a copy of the graph as an image file. An alternative copying option is to click the icon in the top right corner of the chart and drag it into a Microsoft Word or PowerPoint document.

Spreadsheet Options

- Gene expression analysis results are displayed in the spreadsheet (Figure 5).
- Click on any column header to display the data sorted by the column values.
- Right click on the spreadsheet to export the results to a Microsoft Excel spreadsheet or an alternative format.

Excel, Microsoft, and PowerPoint are trademarks of Microsoft Corporation.

Bio-Rad's real-time thermal cyclers are licensed real-time thermal cyclers under Applera's United States Patent Number 6,814,934 B1 for use in research, human in vitro diagnostics, and all other fields except veterinary diagnostics.

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.

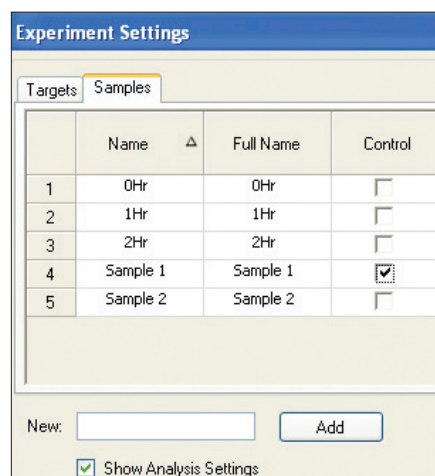


Fig. 3. Assigning the control sample.

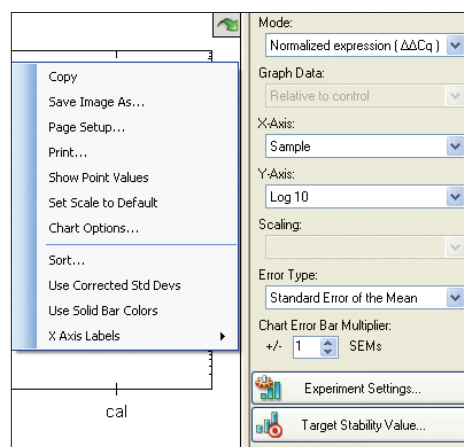


Fig. 4. Graphing analysis options and right-click graph menu.

Target	Sample	Ctl	Expression	Expression SD	Corrected Expression SD
NGK	0Hr	*	1.00000	0.12920	0.12920
NGK	1Hr		2.57905	0.50071	0.50071
NGK	2Hr		18.46978	3.94070	3.94070
NGK	3Hr		163.08526	27.34259	27.34259
NGK	4Hr		1840.02543	250.03713	250.03713
NGK	5Hr		18063.89405	2972.66882	2972.66882

Fig. 5. Gene expression results in spreadsheet format.



**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 31 3689 6600 Canada 905 364 3435 China 86 20 8732 2339 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 30 210 777 4396 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 0508 805 500 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 061 717 95 55 Taiwan 886 2 2578 7189 United Kingdom 020 8328 2000