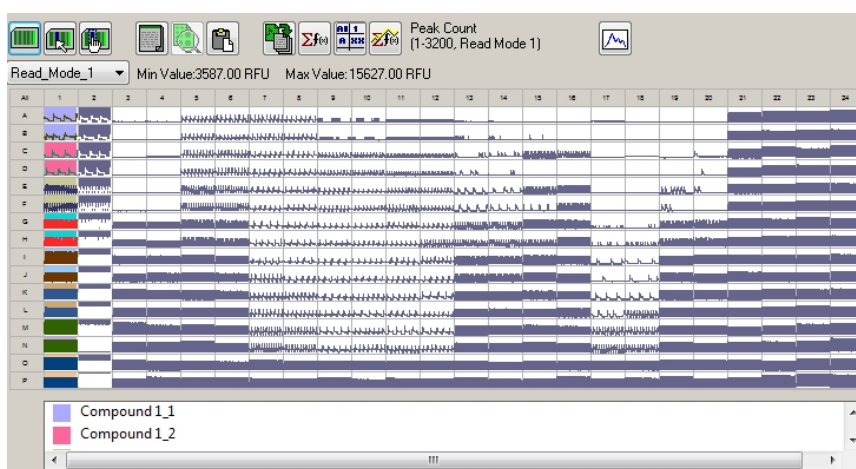



ScreenWorks Peak Pro Software Version 2.0 Quick Start


The ScreenWorks® Peak Pro™ Software Version 2.0 requires the purchase of a separate module license after a 14-day trial. After you have saved experiment data files, you can run the Peak Pro 2.0 software module. The Peak Pro 2.0 software module functionality is best suited for cardiomyocyte, neuronal, and similarly oscillating assay data. To help you understand a little about what this functionality can do for you, this quick start guide uses the provided example cardiomyocyte assay data file called *Demo_Peak_Pro_2.fmd*.

To run the Peak Pro 2.0 software module:

1. Select **File > Open > Data File(s)**.
2. In the **Open** dialog, select your data file name, then click **Open**.
3. In the Multi-Well Graph, select the individual wells, rows, or columns to analyze.



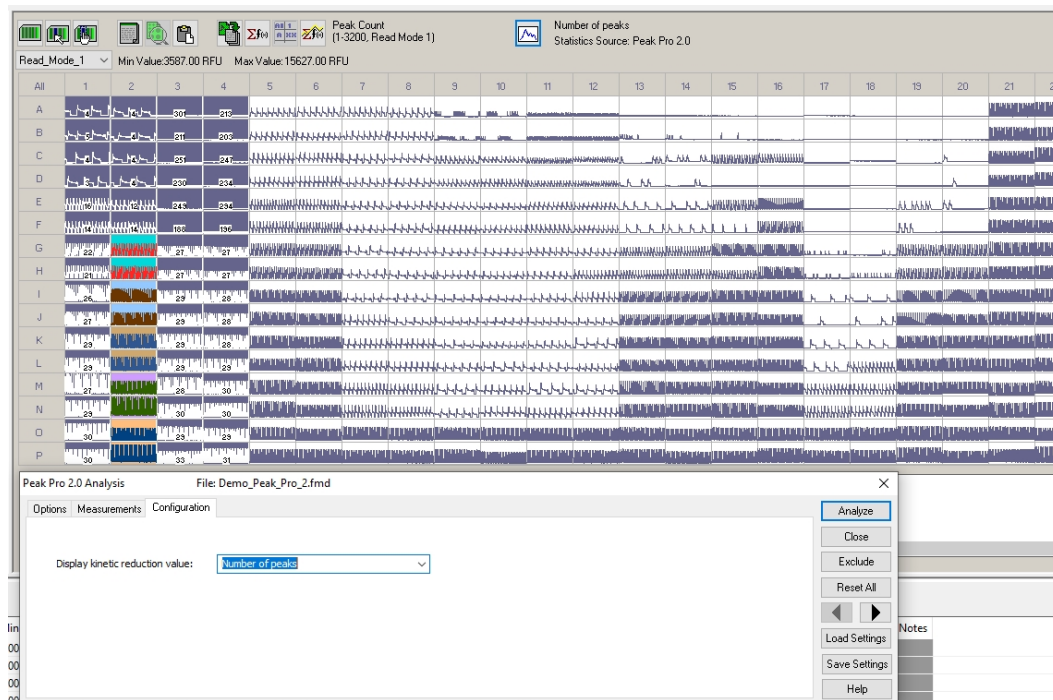
4. Click  to open the **Peak Pro 2.0 Analysis** dialog and collate the selected data, compute automatic baselines, thresholds, optimal search vector length, and format the results for subsequent analysis, sorting and recording.

 **Tip:** If needed, you can change your well selections.

5. In the **Options** tab, either run the default settings or manually adjust.
6. In the **Measurements** tab, if needed, adjust ranges.

 **Note:** All the listed measurements generate during the analysis.

- (Optional) In the **Configurations** tab, select one statistic type to **Display kinetic reduction value** in the Multi-Well Graph for your selected wells, for example, display **Number of peaks**.



- To start the analysis, click **Analyze**.

9. Review the results in the **Detail Graph** pane.

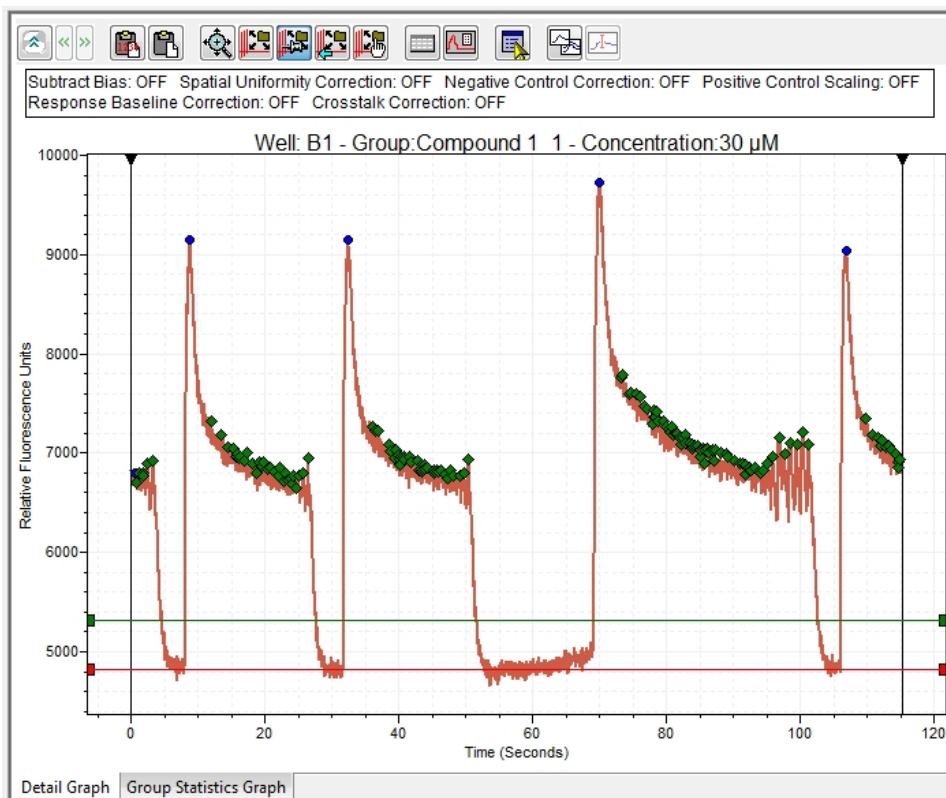
Each major peak is marked with a blue dot. The green diamond markers are secondary peaks.



Tip: When there are a lot displayed, the secondary markers are questionably significant. Adjust the settings so that there is only one major peak and few secondary peaks per event.

The major peak is the first peak in an event above the trigger level.

An event is that portion of a signal that crosses the trigger level in both directions.

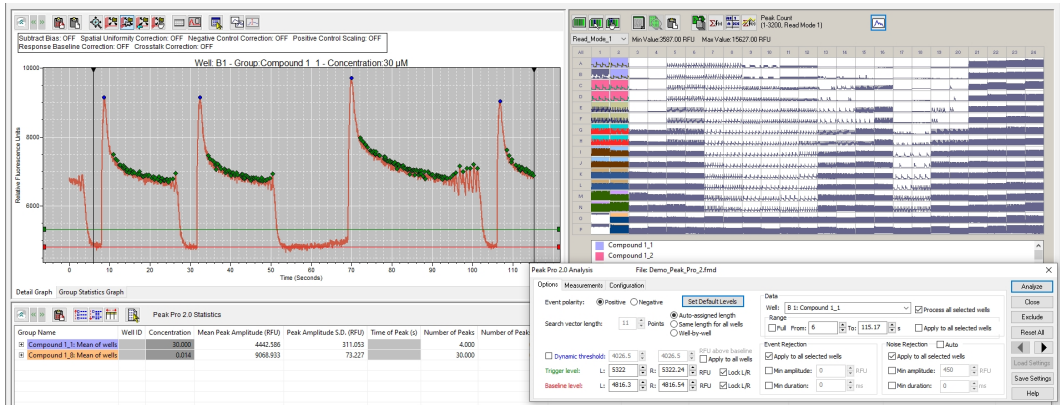


Using the **Peak Pro 2.0 software module** dialog arrow keys, you can scroll through each well graph to examine the result individually. Alternatively, you can select wells directly in the **Options** tab, in the **Data > Well** list.

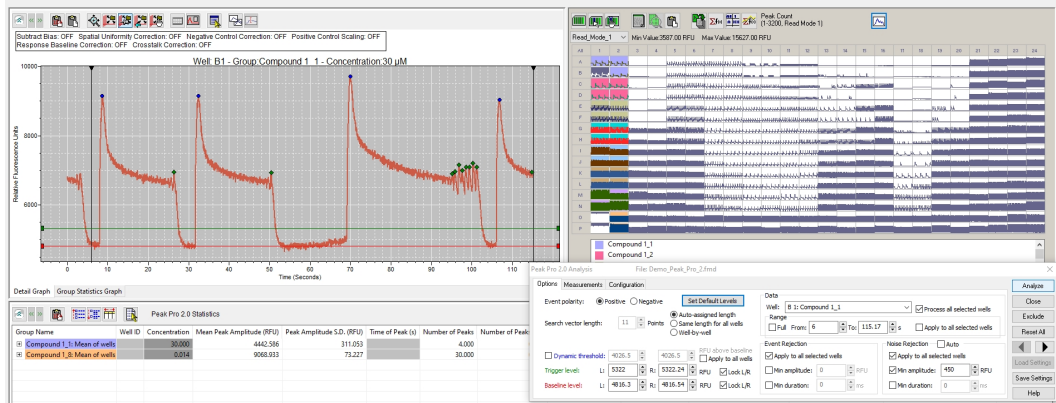
- If the first event marked is a partial signal that you want to skip in your analysis, adjust the analysis data time range.

To adjust the analysis data time range:

- Do one of the following:
 - In the graph, you can manually move the vertical start time range line.
 - In the **Options** tab, in the **Data** section, deselect **Full**, and type a **From** starting value.
- In the **Options** tab, click **Analyze**.




- To reduce the number of secondary peaks, which are likely noise and not biologically significant, adjust settings in the **Options** tab, click **Analyze**, and review the results again in the **Detail Graph** pane.



- After you reduce all of the possible noise and have clearly identified the real biological events, you are finished with the analysis. You can click **Save Settings** to store them with your data file to load later, or just close the floating dialog.



Tip: To reopen the **Peak Pro 2.0 Analysis** dialog, click . Your last settings are retained in the dialog as long as your data file stays open.

Saved settings load automatically the next time the saved data file opens in the Peak Pro 2.0 software module dialog. If you change the settings and want to return to the originally saved settings, click **Load Settings**.

- View the results in the **Peak Pro 2.0 Statistics** pane.

Group Name	Well ID	Number of Peaks	Number of Peaks S.D.	Mean EAD-like peak Rate (PpM)	EAD-like peak Rate S.D. (PpM)	CTD @ 10% (s)	CTD @ 10% S.D.
Compound 1.1: Mean of wells		4.000	0.000			0.702	0.060
Compound 1.8: Mean of wells		29.000	0.000			0.544	0.029

In the statistics pane, the results are sorted by **Group Name**.



Note: **Group Statistics Graph** functionality is not supported in ScreenWorks Software version 5.1.

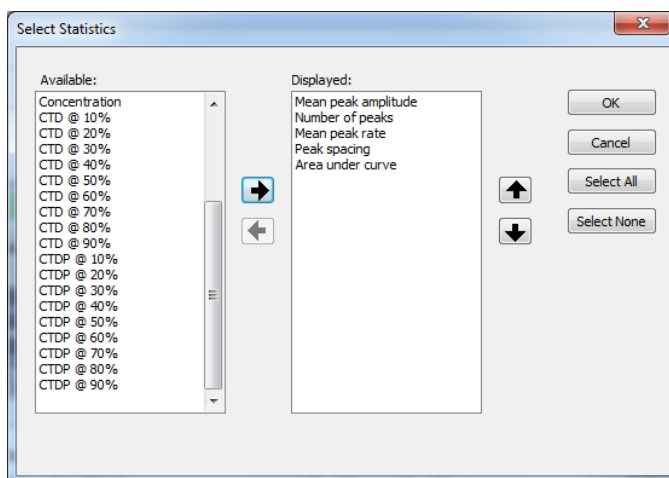
- Expand the **Group** rows and the **Mean** rows to see the **Event** rows of data.

Group Name	Well ID	Number of Peaks	Number of Peaks S.D.	Mean EAD-like peak Rate (PpM)	EAD-like peak Rate S.D. (PpM)	CTD @ 10% (s)	CTD @ 10% S.D.
Compound 1.1: Mean of wells		4.000	0.000			0.702	0.060
B 1: Mean of events	B 1	4				0.702	0.060
Event 1	B 1			170.795	118.585	0.660	
Event 2	B 1			301.752	330.202	0.744	
Event 3	B 1			244.533	249.135		
Event 4	B 1			175.054	148.435		
Compound 1.8: Mean of wells		29.000	0.000			0.544	0.029
O 2: Mean of events	O 2	29				0.544	0.029
Event 1	O 2					0.576	
Event 2	O 2					0.563	
Event 3	O 2					0.537	
Event 4	O 2					0.584	
Event 5	O 2					0.579	
Event 6	O 2					0.563	
Event 7	O 2					0.568	

Each main row contains the average of measurements of each event as selected, and each event row contains the individual metrics of the individual peaks and related measurements.



- To specify which of the data results display in the statistics pane, click **Select Statistics**.
- In the **Select Statistics** dialog, choose which statistics display and in which order.



17. When finished making changes, click **OK**.



Note: The **Peak Pro 2.0 Statistics** pane data never saves with the open data file used for analysis.

18. To save your analysis data, do either of the following:

- In the Multi-Well Graph pane, click  **Export**.
- In the Statistics pane, click  **Copy Table Data to Clipboard** and then paste the delimited data into third-party software of your choice.

Externally, the saved data can be converted into graphs and reports.

Contact Us

Phone: [+1-800-635-5577](tel:+1-800-635-5577)
Web: moleculardevices.com
Email: info@moldev.com

Visit our website for a current listing of worldwide distributors.