

Setup for Z'-LYTE® assay on SpectraMax® Paradigm® Microplate Detection Platform with SoftMax® Pro 6 software

The Molecular Devices SpectraMax® Paradigm® Microplate Detection Platform was tested for compatibility with Life Technologies' Z'-LYTE® FRET assays using the Z'-LYTE® Tyr6 kit (PV4122) against JAK2 JH1/JH2 kinase. The following document is intended to demonstrate setup of this instrument.

For more detailed information and technical support of Life Technologies assays, please call 1-800-955-6288 and enter extension 40266 or email drugdiscoverytech@lifetech.com.

For more detailed information and technical support of Molecular Devices instruments or software, please contact Molecular Devices at 1-800-635-5577 or www.moleculardevices.com.

Setup Guide on the Molecular Devices SpectraMax® Paradigm® Microplate Detection

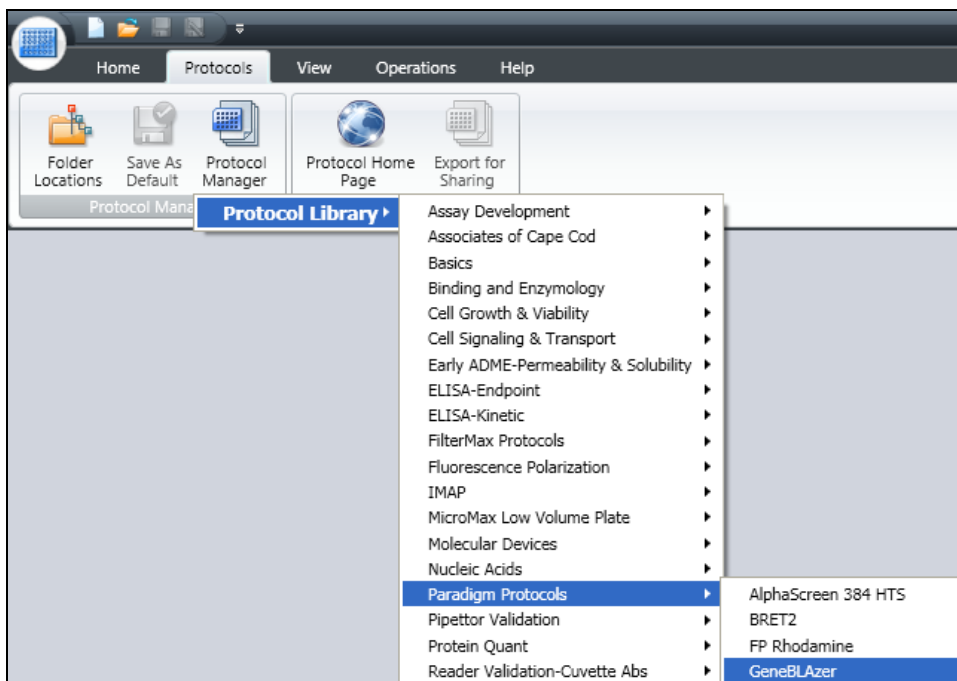
A. Recommended Optics

Parameter	Specification
Detection Cartridge Name	SpectraMax® Paradigm® Fluorescence Intensity (FI) GeneBLAzer® Detection Cartridge
Part Number	0200-7006
Detection Technique	FRET, Fluorescence Intensity
Light Source	LED, ultra high power
Filter Set	EX: 406/15 EM1: 465/35 EM2: 535/25
Applications	Designed for use with GeneBLAzer® reagents; also suitable for Z'-LYTE® assays

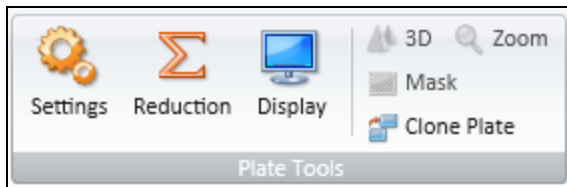
Setup Guide on the Molecular Devices SpectraMax® Paradigm® Microplate Detection

B. Instrument Setup:

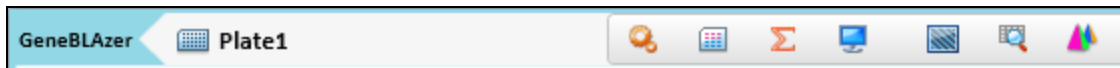
1. Open SoftMax® Pro 6 software. Click on "Protocol Manager" to open the Protocol Library. Within the "Paradigm® Protocols" folder, locate the "GeneBLAzer®" protocol and click to open.



2. Click on "Plate01" in the Navigation Tree on the left side of the screen. Click on the Settings icon either in the toolbar at the top of the screen...

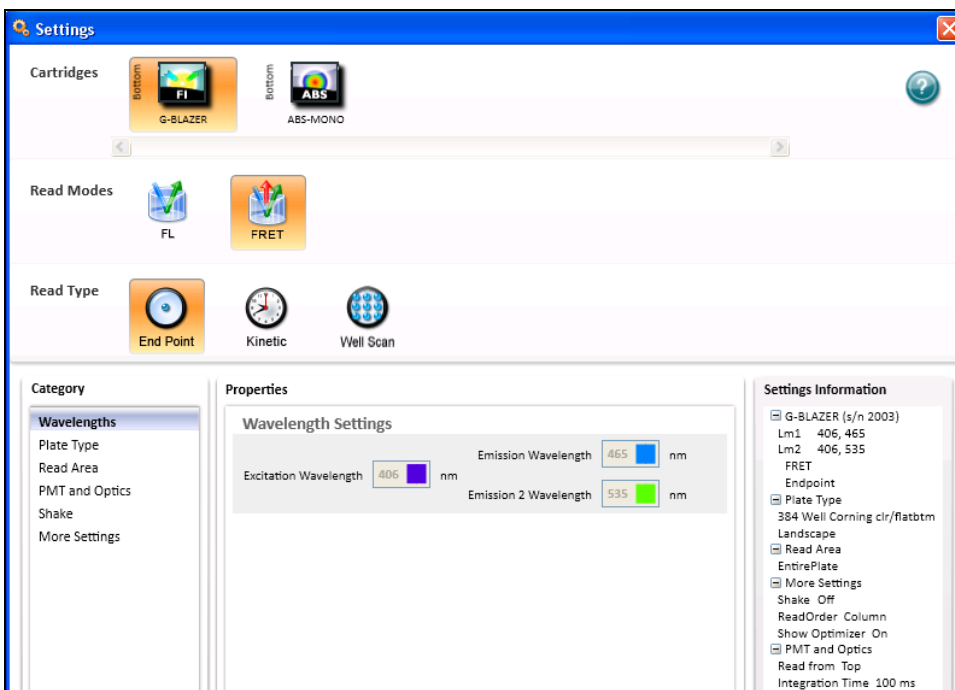


...or in the plate section header.

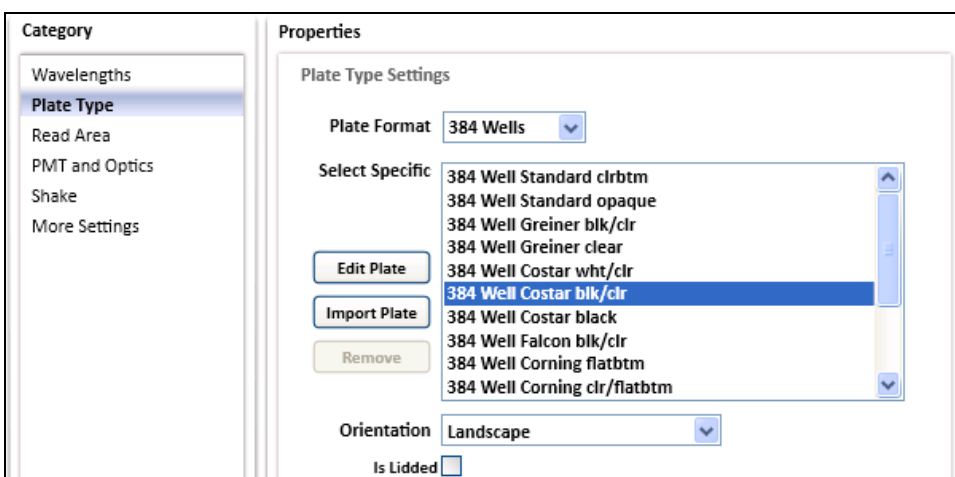


Setup Guide on the Molecular Devices SpectraMax® Paradigm® Microplate Detection

3. This opens the Settings window. The GeneBLAzer® (G-BLAZER) cartridge and its wavelengths already appear under Wavelength Settings.

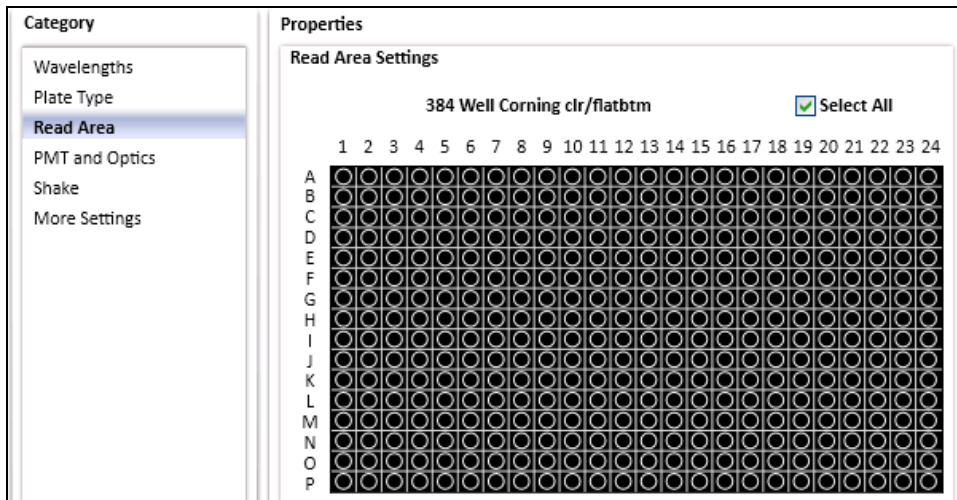


4. Choose the desired plate type, using the upper dropdown menu to choose plate format (96, 384, or 1536 wells) and the "Select Specific" menu to choose the specific plate type.

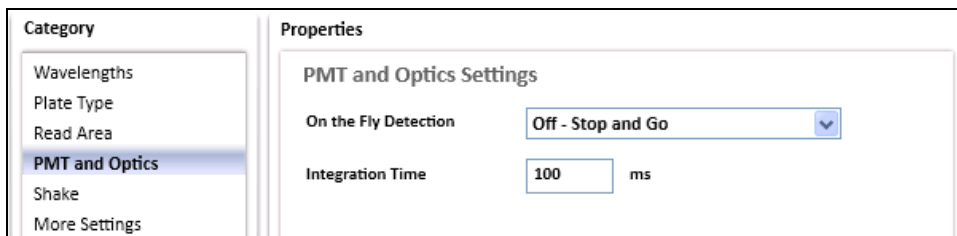


Setup Guide on the Molecular Devices SpectraMax® Paradigm® Microplate Detection

5. Now select the area of the plate to read.



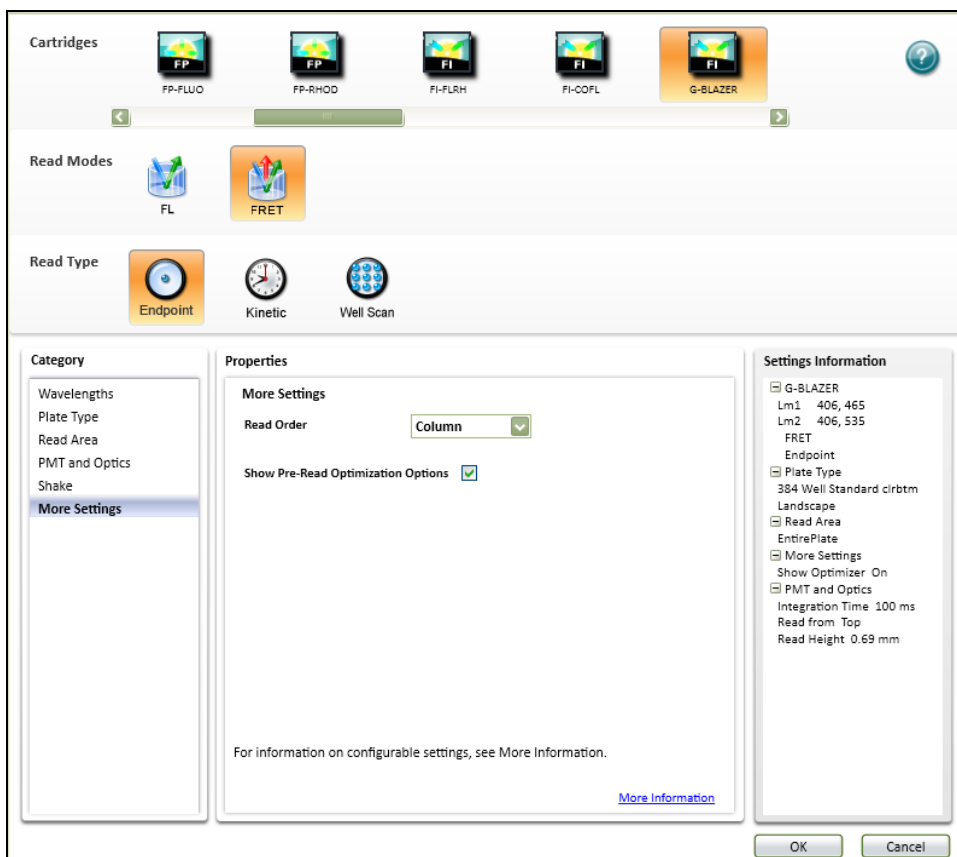
6. PMT and Optics Settings include the option to read using On the Fly detection and adjust the integration time if desired. "Off – Stop and Go" is the default setting. To select On the Fly for faster read times, use the dropdown menu to choose Performance or Speed (faster) On the Fly options. The default integration time is 140 msec. Shorter integration times enable faster reading, while longer integration times enable better performance.



Setup Guide on the Molecular Devices SpectraMax® Paradigm® Microplate Detection

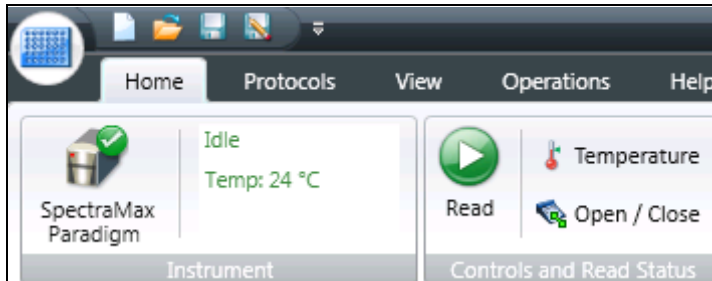
7. In the category "More Settings", choose the read order corresponding to how the assay plate is set up. If the entire plate is to be read, choose "Row". If entire rows of a partial plate are to be read, choose "Row"; if entire columns of a partial plate are to be read, choose "Column". Check the box "Show Pre-Read Optimization Options" to enable the Microplate Optimization and Read Height Adjustment options upon initiation of the plate read.

Click OK to close the Settings window.



Setup Guide on the Molecular Devices SpectraMax® Paradigm® Microplate Detection

8. To read the plate, click the green "Read" button at the top of the screen.

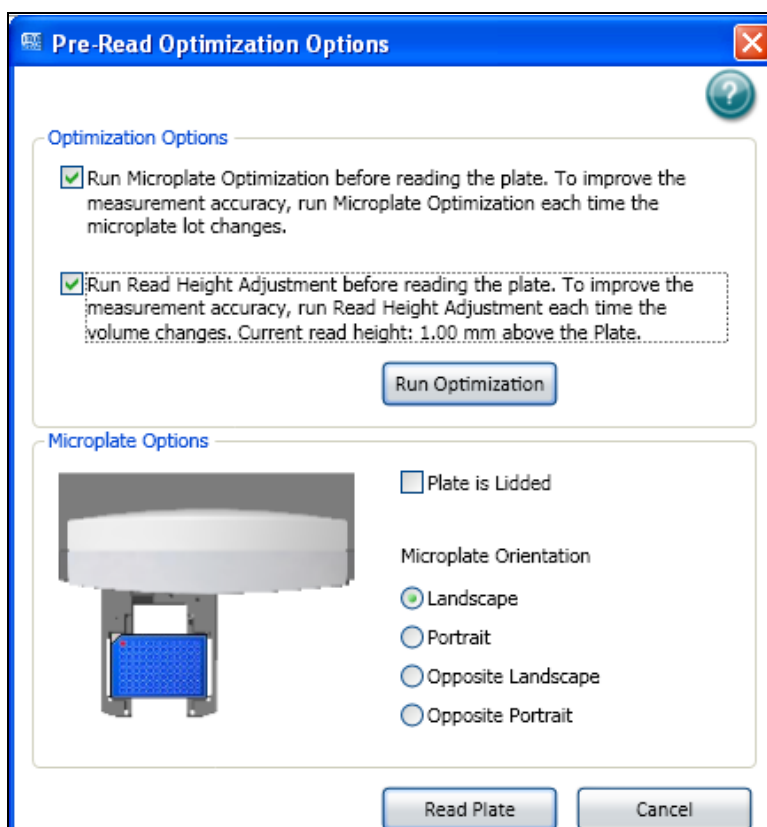


Setup Guide on the Molecular Devices SpectraMax® Paradigm® Microplate Detection

9. If selected, pre-read optimization options will appear:

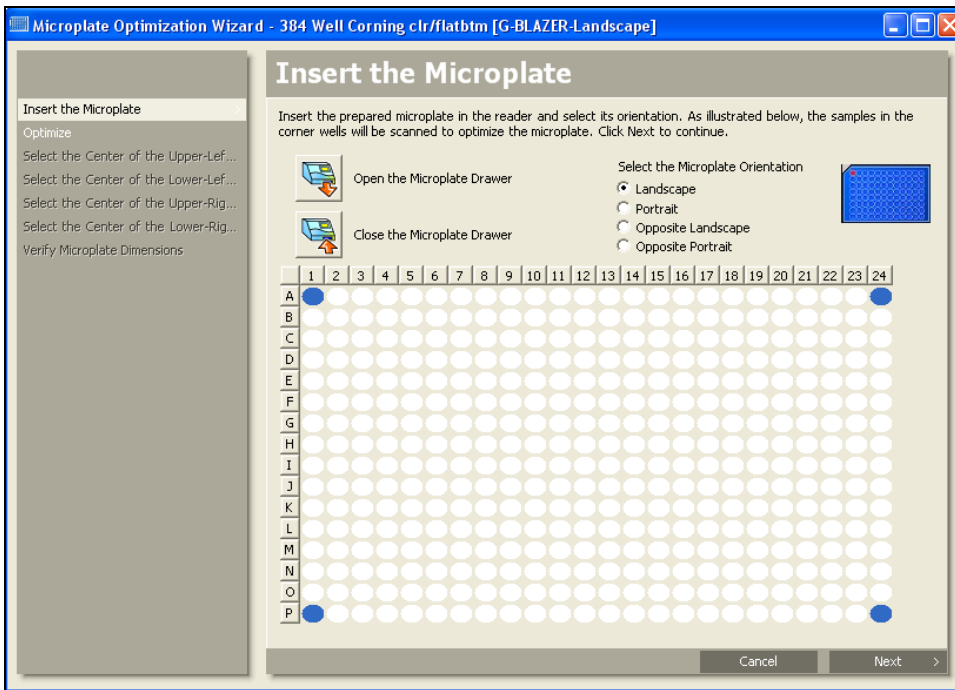
- Microplate Optimization scans the four corner wells of the plate and adjusts the microplate dimensions if necessary to improve accuracy. It requires that all four corners of the microplate contain detectable fluorescent material (i.e. positive control samples).
- Read Height Adjustment determines the height above the plate at which the best signal is detected. It can be performed using any well in the plate with a relatively strong fluorescent signal (i.e. positive control sample).
- If the plate is lidded, check the box. Make sure that the selected microplate orientation matches the orientation of the actual assay plate.

Click "Run Optimization" to proceed. Alternatively, if no optimization is desired, leave the boxes unchecked and click "Read Plate."

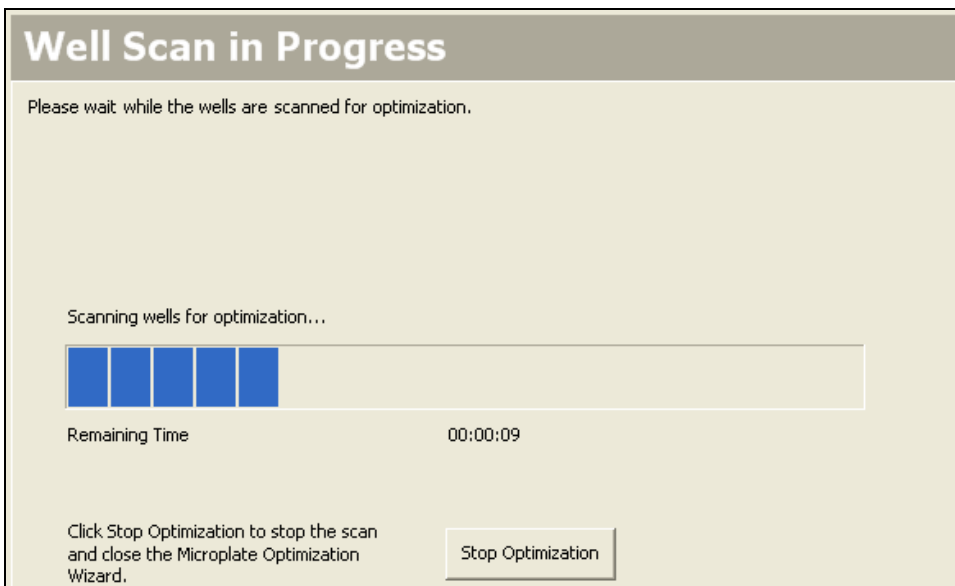


Setup Guide on the Molecular Devices SpectraMax® Paradigm® Microplate Detection

10. If optimization was selected, a wizard will pop up. Follow the steps outlined in the wizard.

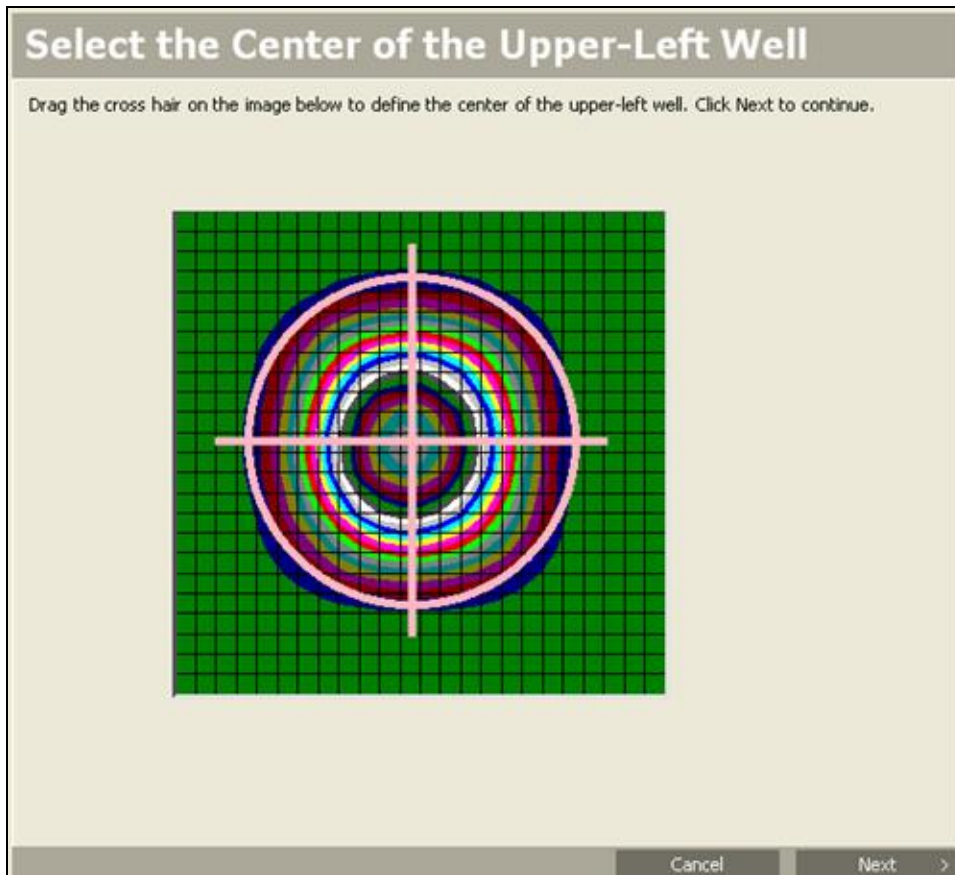


11. When you select read plate, a progress screen will appear as the plate is read.



Setup Guide on the Molecular Devices SpectraMax® Paradigm® Microplate Detection

12. Center the pink target over the image of the scanned well. Click "Next" and repeat for the remaining three wells. This adjusts the microplate definition to match the actual plate.



Setup Guide on the Molecular Devices SpectraMax® Paradigm® Microplate Detection

13. Click "Save" to save the modified plate dimensions with the Microplate Name as shown. This optimized microplate type will be available in the Settings for future use.

Verify Microplate Dimensions

Verify the dimensions of the microplate. You can edit the values in the fields or return to a well step to redefine its center. Type a name for the microplate definition in the Microplate Name field. Click Save to save the microplate definition.

<input type="checkbox"/> Microplate Dimensions	
Bottom-row y offset (mm)	8.99
Column spacing (mm)	4.5
Left-column x offset (mm)	12.12
Right-column x offset (mm)	12.12
Row spacing (mm)	4.5
Top-row y offset (mm)	8.99

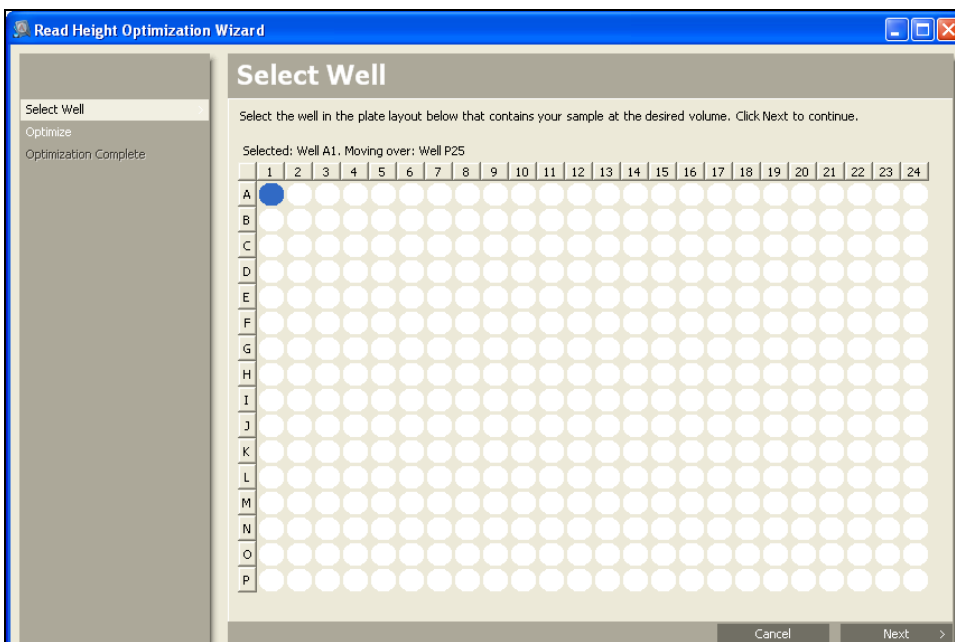
<input type="checkbox"/> Microplate Name	
Microplate Name	384 Well Corning clr/flatbtm [G-BLAZER-Landscape]

Bottom-row y offset (mm)
The distance in millimeters from the lower edge of the microplate to the horizontal center of the bottom row.

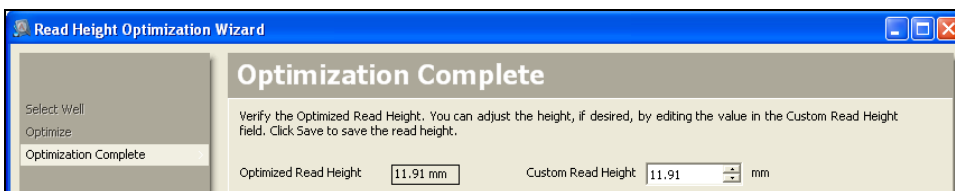
Cancel
< Back
Save

Setup Guide on the Molecular Devices SpectraMax® Paradigm® Microplate Detection

14. If you chose to perform Read Height Adjustment, this wizard will now appear. Select the well you want to use for read height adjustment. This should be a relatively bright well, e.g. a positive control. Click "Next" to read.

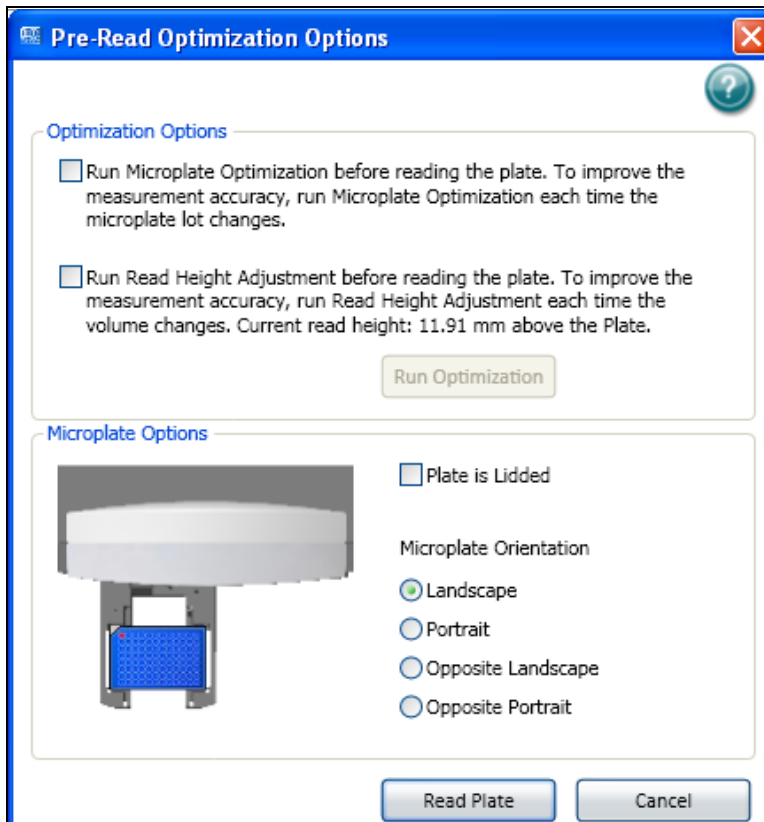


15. The instrument will calculate and report optimized read height.



Setup Guide on the Molecular Devices SpectraMax® Paradigm® Microplate Detection

16. After optimization is complete, click on "Read Plate" to proceed.



Setup Guide on the Molecular Devices SpectraMax® Paradigm® Microplate Detection

17. After the plate is read, data will appear in the plate section:

The screenshot shows the SpectraMax software interface with the following components:

- Navigation Tree:** Shows a hierarchy starting with 'Expt1', followed by 'Plate1', and then various assay parameters like 'JAK2 JH1/JH2', 'V617F', '0%phos', '100%phos', 'JAK2_inh', 'V617F_inh', 'Background', and 'Graph1'.
- Document/Comparison Tab:** The main window title is 'Expt1 Plate1'. A toolbar contains icons for search, print, sum, and other functions.
- Plate1 Data Table:** A grid with columns 1-24 and rows A-P. The data values are:

	1	2	3	4	5	6	7	8	9	10	11	12	13	14	15	16	17	18	19	20	21	22	23	24
A	4e7	2e5	9e7	9e7	9e7	9e7	9e7	9e7	9e7	9e7	9e7	9e7	9e7	9e7	9e7	9e7	9e7	9e7	9e7	9e7	9e7	9e7	9e7	9e7
B	3e7	1e5	2e7	1e7	1e7	9e6	9e6	1e7	1e7	1e7	6e4	1e7	1e7	1e7	9e6	1e7	1e7	1e7	1e7	1e7	7e5	3e6	7e7	9e6
C	3e7	6e4	2e7	1e7	1e7	1e7	1e7	1e7	1e7	1e7	7e4	3e4	2e7	7e6	1e7	1e7	1e7	1e7	1e7	1e7	1e7	4e5	8e6	9e6
D	4e7	2e5	9e7	9e7	9e7	9e7	9e7	9e7	9e7	9e7	1e8	9e5	2e5	6e7	8e7	9e7	9e7	9e7	9e7	9e7	9e7	9e7	9e7	9e7
E	3e7	8e4	3e7	1e7	1e7	1e7	1e7	1e7	1e7	1e7	1e5	3e4	2e7	9e6	9e6	9e6	9e6	9e6	9e6	9e6	9e6	9e6	9e6	9e6
F	6e7	3e5	5e7	8e7	8e7	8e7	8e7	8e7	8e7	8e7	1e6	2e5	7e7	9e7	1e8	1e8	1e8	1e8	1e8	1e8	1e8	9e4	1e8	
G	4e7	2e5	2e7	1e7	9e6	1e7	9e6	1e7	9e6	1e7	2e5	2e4	3e7	1e7	1e7	1e7	1e7	1e7	1e7	1e7	1e7	1e7	2e4	1e7
H	5e7	3e5	6e7	9e7	9e7	9e7	9e7	9e7	9e7	9e7	1e6	3e5	6e7	8e7	9e7	9e7	9e7	9e7	9e7	9e7	9e7	9e7	9e7	9e4
I	4e7	1e5	3e7	1e7	1e7	1e7	1e7	1e7	1e7	1e7	1e5	3e4	2e7	1e7	1e7	1e7	1e7	1e7	1e7	1e7	1e7	1e7	1e4	1e7
J	5e7	3e5	6e7	9e7	9e7	9e7	9e7	9e7	9e7	9e7	1e6	3e5	6e7	8e7	9e7	9e7	9e7	9e7	9e7	9e7	9e7	9e7	9e7	9e4
K	4e7	1e5	2e7	1e7	1e7	1e7	1e7	1e7	1e7	1e7	1e5	3e4	2e7	1e7	1e7	1e7	1e7	1e7	1e7	1e7	1e7	1e7	1e4	1e7
L	4e7	2e5	9e7	9e7	9e7	9e7	9e7	9e7	9e7	9e7	1e8	9e5	2e5	6e7	8e7	9e7	9e7	9e7	9e7	9e7	9e7	9e7	9e7	9e7
M	3e7	9e4	3e7	1e7	2e7	2e7	1e7	1e7	1e7	1e7	2e5	3e4	3e7	1e7	1e7	1e7	1e7	1e7	1e7	1e7	1e7	1e7	1e4	8e6
N	4e7	2e5	9e7	9e7	9e7	9e7	9e7	9e7	9e7	9e7	1e6	2e5	6e7	8e7	9e7	9e7	9e7	9e7	9e7	9e7	9e7	9e7	9e7	9e4
O	4e7	2e5	9e7	9e7	9e7	9e7	9e7	9e7	9e7	9e7	1e6	2e5	6e7	8e7	9e7	9e7	9e7	9e7	9e7	9e7	9e7	9e7	9e7	9e4
P	4e7	1e5	2e7	1e7	1e7	1e7	1e7	1e7	1e7	1e7	1e5	3e4	2e7	1e7	1e7	1e7	1e7	1e7	1e7	1e7	1e7	1e7	1e4	1e7
- Settings Information:** Includes 'G-BLAZER (s/n 2003)', 'Lm1 406, 465', 'Lm2 406, 535', 'FRET', 'Endpoint', 'More Settings', 'Shake Off', 'ReadOrder Column', 'Show Optimizer On', 'PMT and Optics', 'Read from Top', and 'Integration Time 100 ms'.
- Read Information:** Shows 'SpectraMax Paradigm', 'ROM v1.2 b5.1 20.05.2011', 'Start Read : 8:21 AM', '8/12/2011', and 'Mean Temperature : 25 °C'.
- Reduction Settings:** 'Wavelength Combination : 1Lm1'.

18. To set up the template for data analysis, click on Template Editor icon:

The screenshot shows the SpectraMax software interface with the following components:

- Navigation Tree:** Similar to the previous screenshot, showing 'Expt1' and 'Plate1' with various assay parameters.
- Document/Comparison Tab:** The main window title is 'Expt1 Plate1'. The toolbar now has a 'Template Editor' icon highlighted with a yellow box.
- Plate1 Data Table:** The same data table as in the previous screenshot, showing values for columns 1-24 and rows A-P.
- Settings Information:** Similar to the previous screenshot, showing assay parameters and settings.
- Read Information:** Similar to the previous screenshot, showing read details.
- Reduction Settings:** Similar to the previous screenshot, showing 'Wavelength Combination : 1Lm1'.

Have a question?

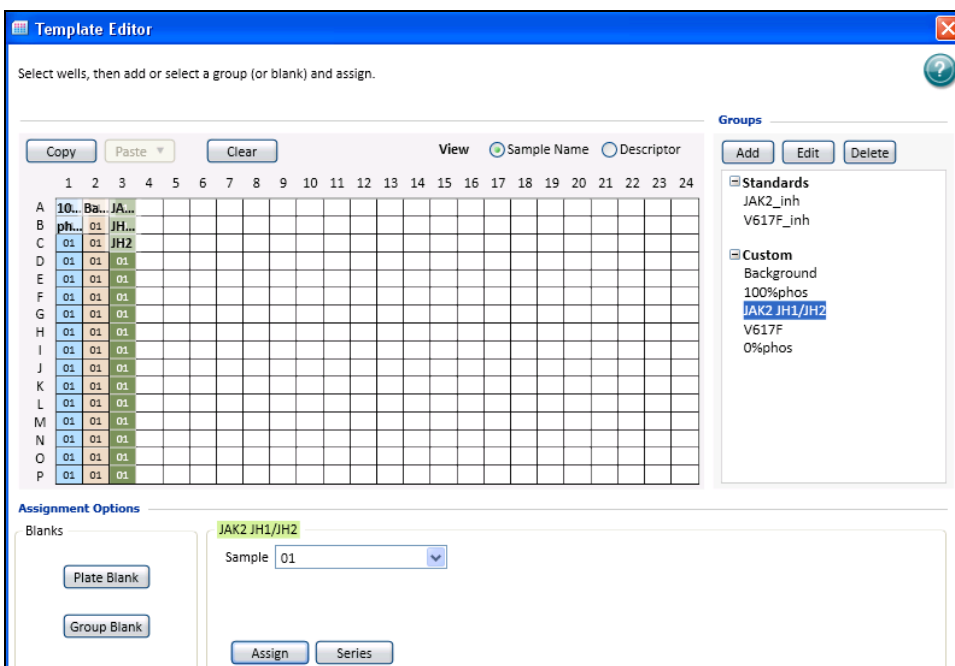
NA: 800-955-6288 or INTL: 760-603-7200 ext. 40266

Contact our Technical Support Team

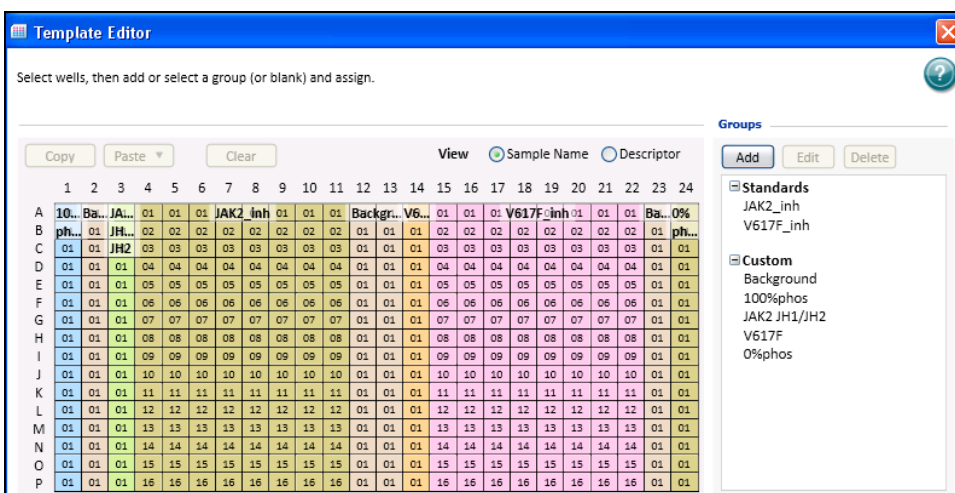
Email: drugdiscoverytech@lifetech.com

Setup Guide on the Molecular Devices SpectraMax® Paradigm® Microplate Detection

19. Select wells and choose the template group you want to assign them to; click Assign. Repeat for each sample type.

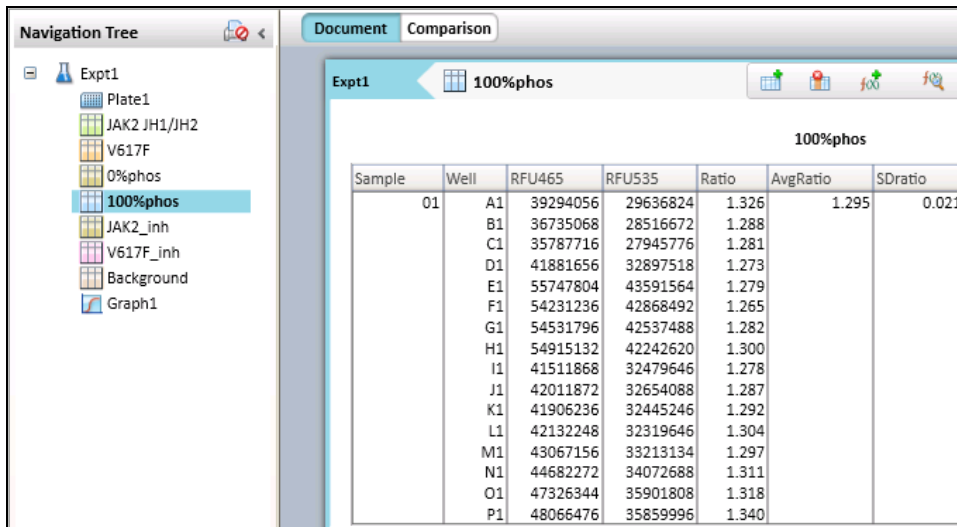


Template with wells assigned to different template groups:



Setup Guide on the Molecular Devices SpectraMax® Paradigm® Microplate Detection

20. When wells are assigned to template groups, data will populate group tables where analysis can be done:



The screenshot shows a software interface with a 'Navigation Tree' on the left and a 'Document Comparison' window on the right. The '100%phos' template group is selected in the tree. The main window displays a data table for this group.

Sample	Well	RFU465	RFU535	Ratio	AvgRatio	SDratio
01	A1	39294056	29636824	1.326	1.295	0.021
	B1	36735068	28516672	1.288		
	C1	35787716	27945776	1.281		
	D1	41881656	32897518	1.273		
	E1	55747804	43591564	1.279		
	F1	54231236	42868492	1.265		
	G1	54531796	42537488	1.282		
	H1	54915132	42242620	1.300		
	I1	41511868	32479646	1.278		
	J1	42011872	32654088	1.287		
	K1	41906236	32445246	1.292		
	L1	42132248	32319646	1.304		
	M1	43067156	33213134	1.297		
	N1	44682272	34072688	1.311		
	O1	47326344	35901808	1.318		
	P1	48066476	35859996	1.340		

Have a question?

NA: 800-955-6288 or INTL: 760-603-7200 ext. 40266

Contact our Technical Support Team

Email: drugdiscoverytech@lifetech.com

Setup Guide on the Molecular Devices SpectraMax® Paradigm® Microplate Detection

C. Results

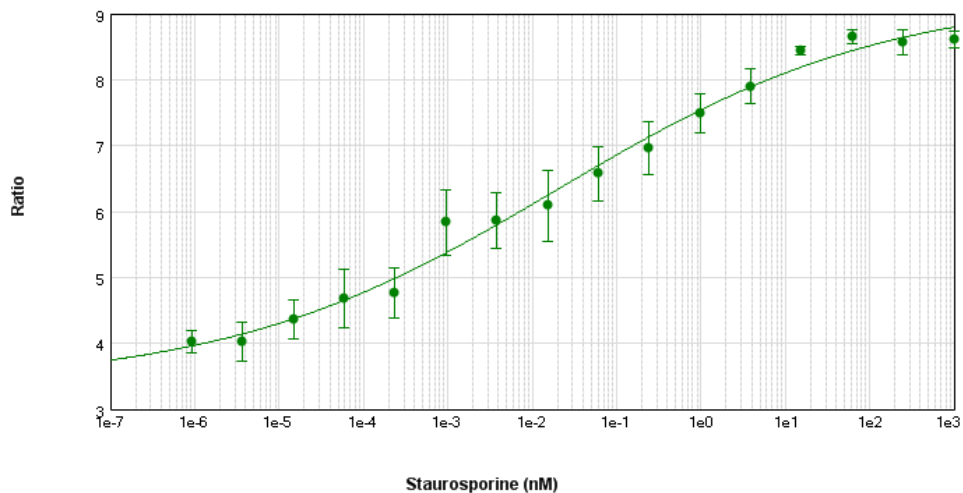


Figure 1: Z'-LYTE® Assay. JAK2 JH1/JH2 Dose-Response Curve read on the Molecular Devices SpectraMax® Paradigm® Microplate Detection Platform. Z' = 0.89.