

LanthaScreen® Terbium Assay Setup Guide on the BioTek Instruments Synergy™ 2 Multi-Mode Microplate Reader

IMPORTANT INFORMATION

Test your plate reader set-up before using LanthaScreen® Terbium and Europium assays

We have developed two technical notes which provide a method for verifying that a fluorescent plate reader is able to detect a change in time-resolved fluorescence energy transfer (TR-FRET) signal, confirming proper instrument set-up and a suitable response. The method is independent of any biological reaction or equilibrium and uses reagents that are on-hand for the LanthaScreen® assay.

For complete instructions, visit www.lifetechnologies.com/instrumentsetup and click on "[Download Terbium assay application note](#)" or "[Download Europium assay application note](#)".

The BioTek Instruments Synergy™ 2 Multi-Mode Microplate Reader was tested for compatibility with Life Technologies LanthaScreen® Terbium-based TR-FRET Assay using the LanthaScreen® fluorescein-Poly GT (PV3610) and Tb-PY20 antibody (PV3552) against JAK2 JH1/JH2 and JAK2 JH1/JH2 V617F kinases. The following document is intended to demonstrate setup of this instrument and provide representative data.

For more detailed information and technical support of Life Technologies assays including specific conditions for assay windows between 2-3 fold, please call 1-800-955-6288 and enter extension 40266 or email drugdiscoverytech@lifetech.com.

For more detailed information and technical support of BioTek instruments or Gen5 software, please call 1-888-451-5171.

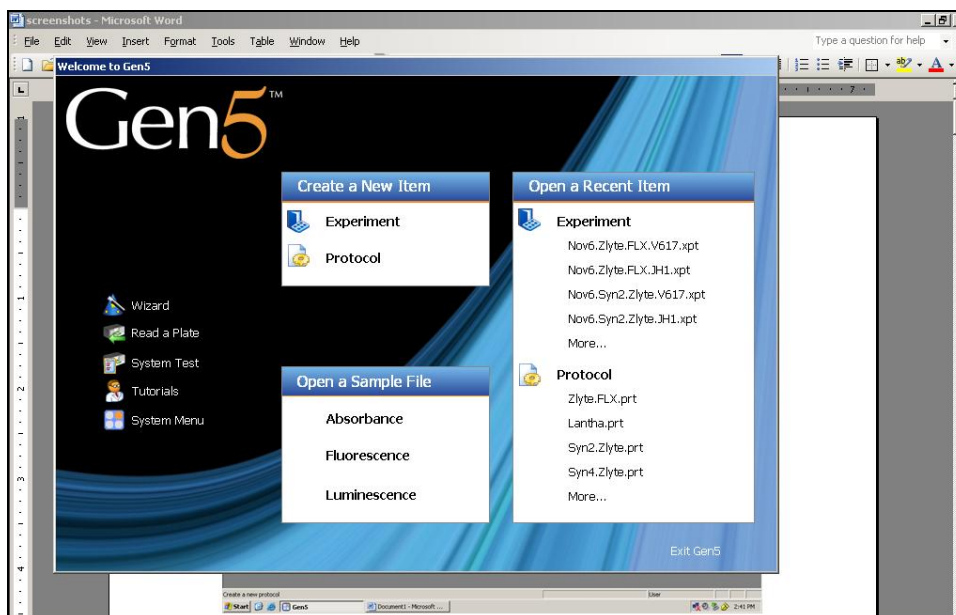
Setup Guide on the BioTek Synergy™ 2 Multi-Mode Microplate Reader

A. Recommended Optics

BioTek Instruments part number	Wavelength (nm)	Diameter (mm)
Excitation (7082230)	340/30	18
Emission 1 (7082268)	495/10	18
Emission 2 (7082267)	520/25	18
Dichroic Mirror (7138400)	400	

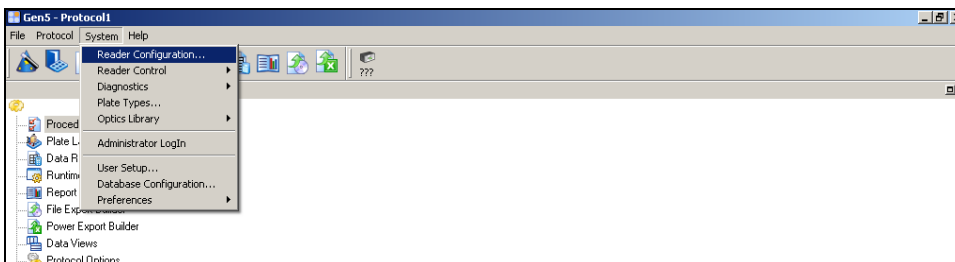
B. Instrument Setup

1. Make certain plate reader is turned on, and open up BioTek Gen5 software on computer.
2. When Gen5 opens, if you do not have a pre-existing protocol for LanthaScreen®, select "Protocol" in the "Create a New Item" menu near the center of the screen.

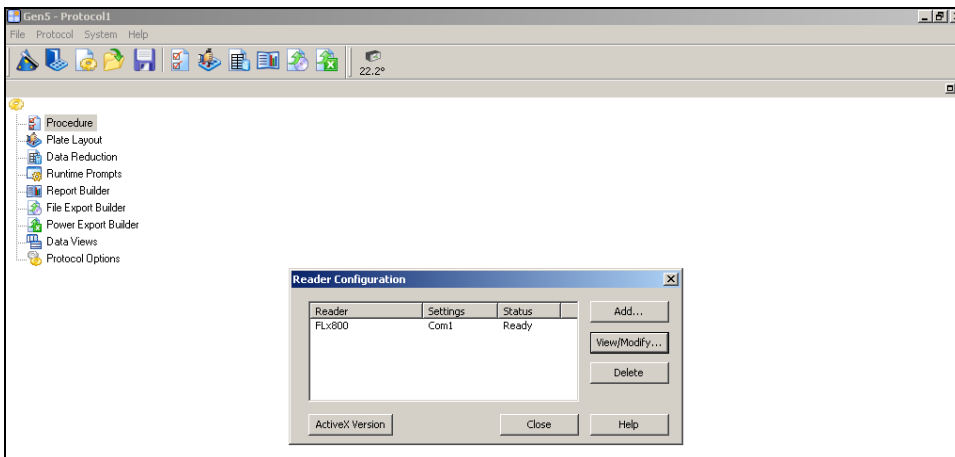


Setup Guide on the BioTek Synergy™ 2 Multi-Mode Microplate Reader

3. At this point, a blank protocol will open (below). To ensure that the correct instrument is selected, click on "Reader Configuration" under tools.

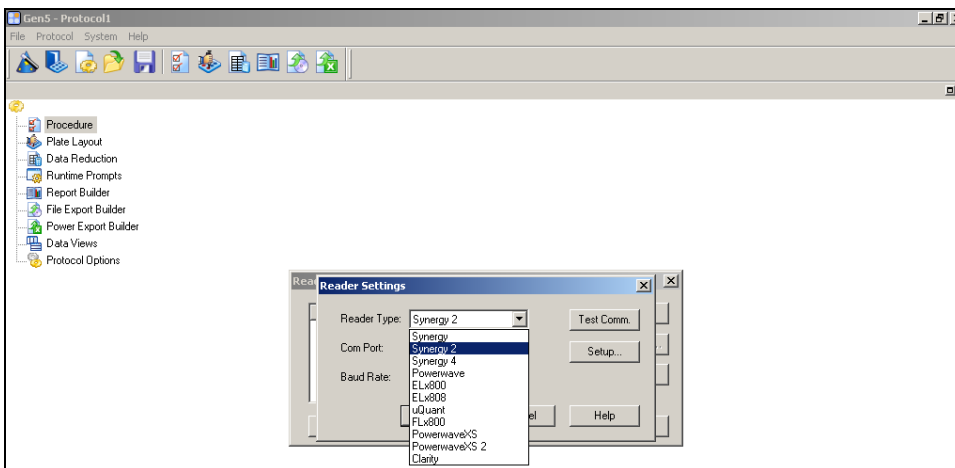


4. Check that the instrument selected is correct (below). If not, select "Delete" to remove the existing instrument, and then "Add".

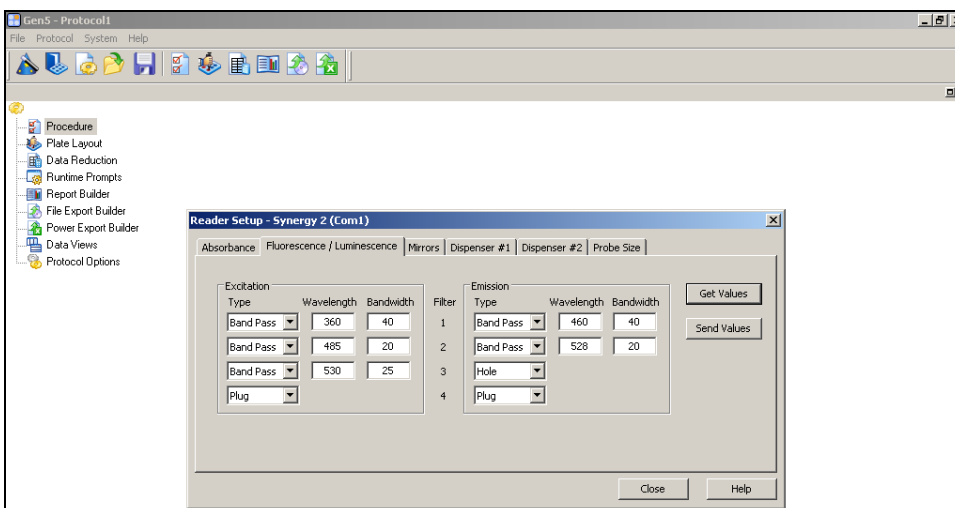


Setup Guide on the BioTek Synergy™ 2 Multi-Mode Microplate Reader

5. Upon selecting "Add" a popup window will appear allowing you to select your instrument from the drop-down menu. Select your instrument here. Click on "Test Communication" to verify that the instrument is communicating properly with the software. When this is done, click on "Setup" finally to check filter and dichroic settings.



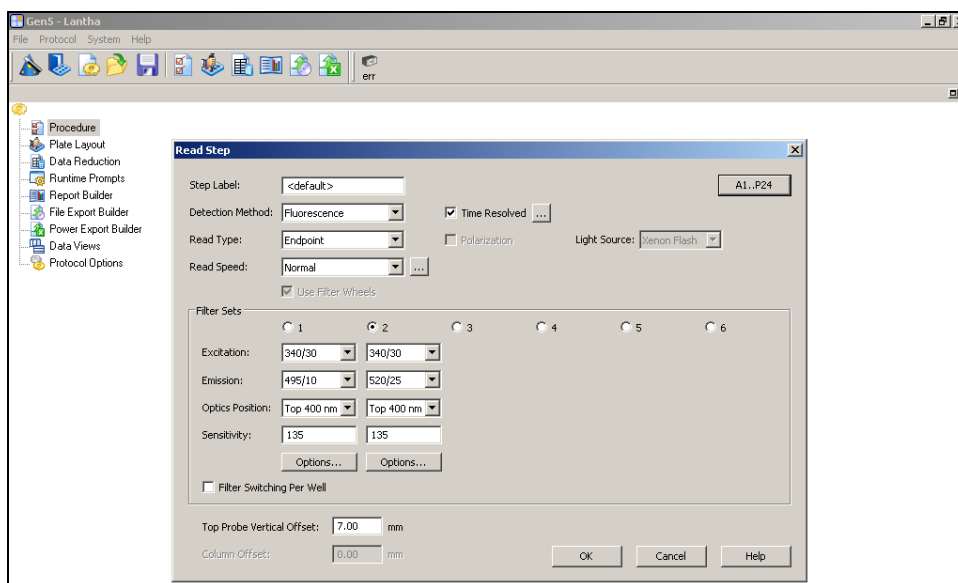
6. Select the "Fluorescence/Luminescence" tab at the top of the popup window. Enter the correct values for your filters present in your filter holder(s). Make certain to do so in the proper positions. Note if you have more than one filter holder for excitation or emission you may have to check that the correct set is listed in the settings or enter the appropriate one. Click on "Send Values" to load the filter positions and settings into the instrument.



Setup Guide on the BioTek Synergy™ 2 Multi-Mode Microplate Reader

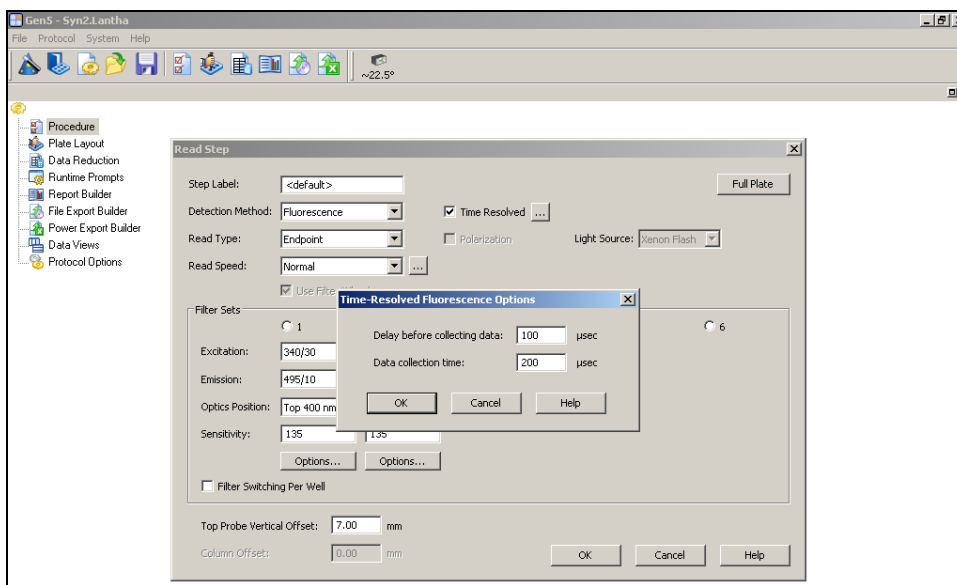
7. Select "Mirrors" and repeat for the dichroic mirror as well, if needed. Select "Close" when finished.

8. Since you have already selected to open a new protocol, at this point the instrument and software are ready to set up a new protocol. Select "Procedure" on the left hand side and a new window will open up. Select "Fluorescence" and "Endpoint" from the drop-down menus. Select the proper filters and dichroic mirror settings, as shown below. **Note:** LanthaScreen® emission filters can be critical, and failure to use appropriate filters can compromise assay results.



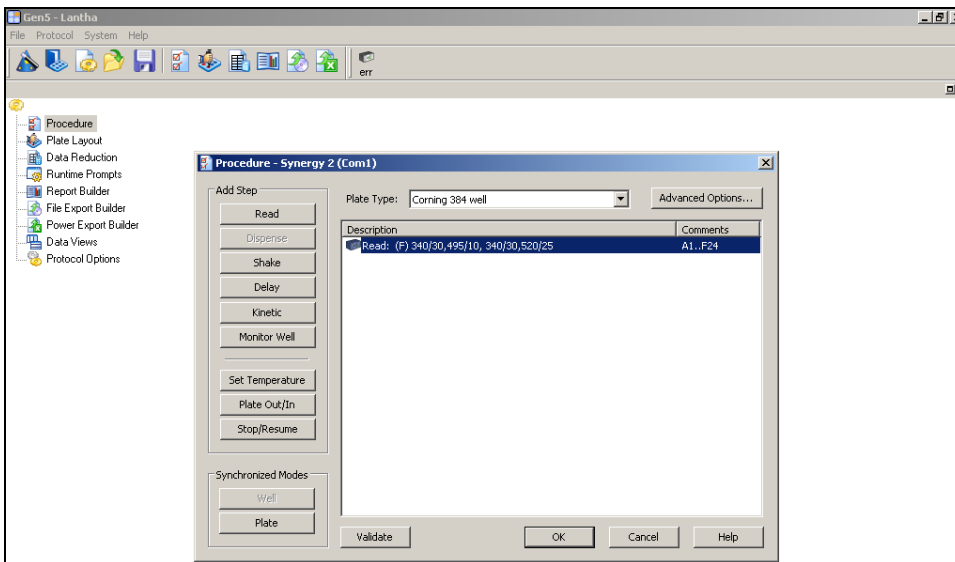
Setup Guide on the BioTek Synergy™ 2 Multi-Mode Microplate Reader

9. Since LanthaScreen® is a Time Resolved FRET assay, select "Time Resolved" in the checkbox to the right of the Detection Method menu. A smaller popup window will appear now; we recommend setting the delay at 100 μ sec and the collection time at 200 μ sec. Click "OK" and then click "OK" again in the main Read Step window.

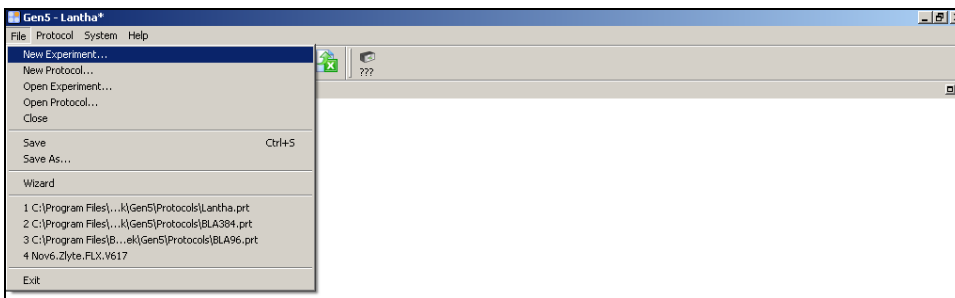


Setup Guide on the BioTek Synergy™ 2 Multi-Mode Microplate Reader

10. If you haven't already, select plate size now. **Note:** we have selected for a Corning 384-well. A "Read" step will appear in blue. Select "Validate" to ensure the protocol steps are valid with the current instrument settings. If the sequence is valid, select "OK" both in the smaller popup window and in the main screen. Select "Save As" from the "File" menu at the top of the screen, and save the protocol with an appropriate name to preserve these instrument settings.

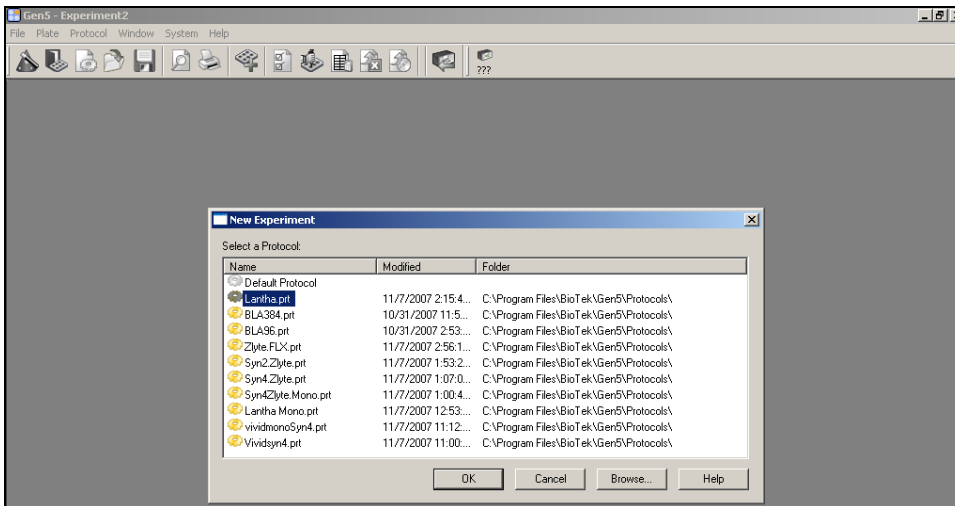


11. To run an actual experiment, click on "File" again, and select "New Experiment" from the drop-down list.

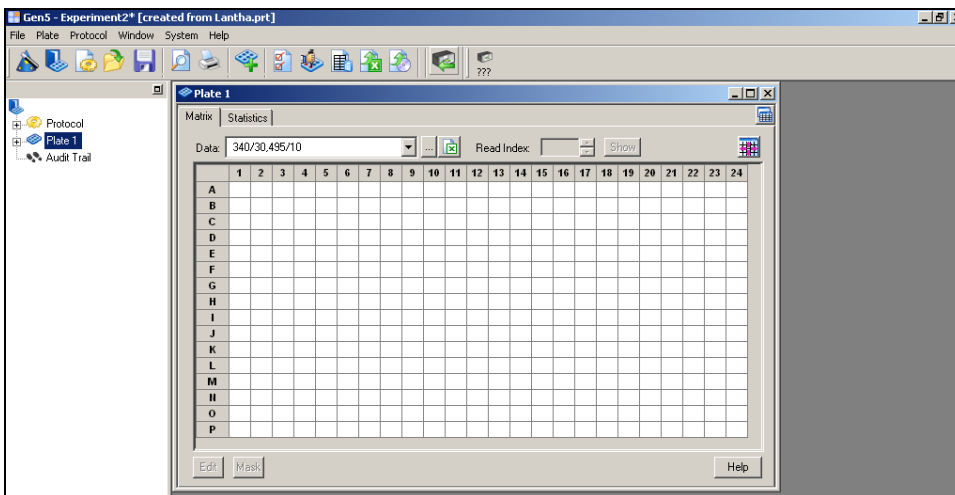


Setup Guide on the BioTek Synergy™ 2 Multi-Mode Microplate Reader

12. A smaller window will appear; select your protocol and click "OK".

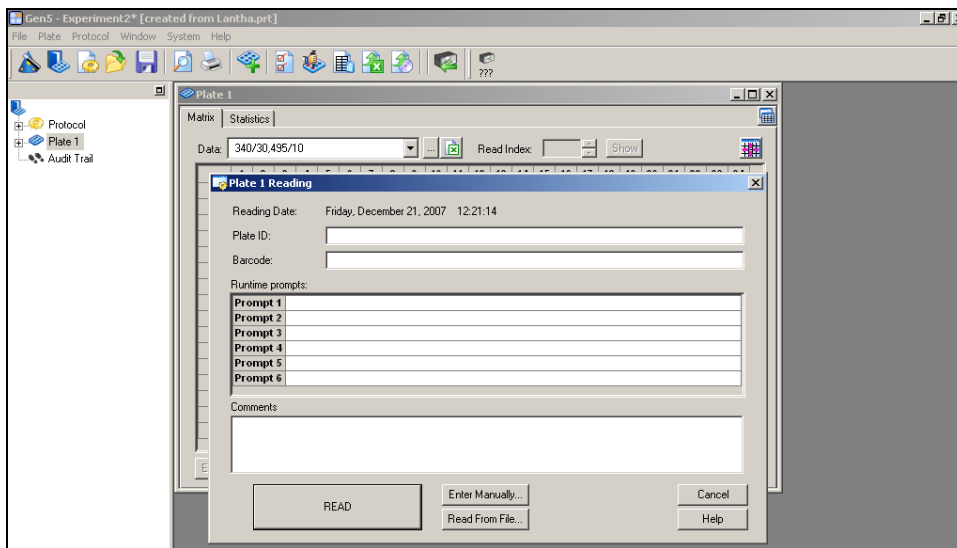


13. A new window will again appear; this is the experiment window. Insert your plate to be read and click on the "Read Plate" icon with the small green arrow (second from last icon in the lower toolbar).



Setup Guide on the BioTek Synergy™ 2 Multi-Mode Microplate Reader

14. One last popup will appear. Select "Read" and instrument will commence reading your plate.
Note: There are several options for collecting data from the BioTek Gen5 software. Data can be exported directly to excel, or appropriate macros can be prepared within the BioTek software. For more information on data output options, please contact BioTek Technical Support.



C. LanthaScreen® Kinase Assay using JAK2 JH1/JH2 and JAK2 JH1/JH2 V617F

Note: The following is a sample titration assay performed for demonstration purposes. The instrument settings above would be sufficient for any LanthaScreen® assay, the information below is provided as representative data. Assay was run at ATP $K_{m_{app}}$ and a kinase concentration producing approximately 70-80% of maximal phosphorylation. ATP and kinase concentrations should be optimized for each kinase by the actual user and titrations/plate layout may be optimized as well. For more information on setting up assays, visit www.lifetechnologies.com/lanthascreen.

1. Prepare initial 100X serial dilution curves in rows A, E, I, and M of a 384-well plate (Figure 1): Dilute Staurosporine, JAK2 Inhibitor 2, JAK3 Inhibitor, and AG-490 to a 100X initial concentration (100uM) in 100% DMSO. Prepare a set of 1:1 serial dilutions from the initial concentration in a 384-well plate, starting with 80 μ l in column 1 and 40 μ l DMSO in wells 2-20. Add 40 μ l from well 1 to well 2, and then mix well 2, and take 40 μ l from well 2 and add to well 3, mix, and so on.

		1	2	3	4	5	6	7	8	9	10	11	12	13	14	15	16	17	18	19	20	21	22	23	24
Staur.	100X	A	█	█	█	█	█	█	█	█	█	█	█	█	█	█	█	█	█	█	█	█	█	█	█
	4X	B	█	█	█	█	█	█	█	█	█	█	█	█	█	█	█	█	█	█	█	█	█	█	█
		C																							
		D																							
JAK2 Inh. II	100X	E	█	█	█	█	█	█	█	█	█	█	█	█	█	█	█	█	█	█	█	█	█	█	█
	4X	F	█	█	█	█	█	█	█	█	█	█	█	█	█	█	█	█	█	█	█	█	█	█	█
		G																							
		H																							
	I																								
	J																								
	K																								
	L																								
	M																								
	N																								
	O																								
	P																								

Figure 1: Schematic of initial compound dilution. Staurosporine, JAK2 Inhibitor II, JAK3 Inhibitor, and AG-490 were all titrated from a 100 μ M starting concentration in the initial dilution series by preparing a 1:1 dilution curve in DMSO. A secondary dilution to 4X was then prepared in the rows below the initial dilution curve (lighter gray) using kinase buffer.

2. The 100X serial dilution set is then diluted to a 4X working concentration in Kinase Buffer (PV3189, 50 mM HEPES pH 7.5, 0.01% BRIJ-35, 10 mM MgCl₂, 1 mM EGTA) in the row below by adding 2 μ l of diluted inhibitor from the well above to 48 μ l of kinase buffer. This will produce a final serial dilution starting at 4 μ M, which will then produce a final assay concentration starting at 1 μ M.

3. Begin to prepare an assay plate: Add 2.5 μ l of the compound dilutions per well into a low volume NBS, 384-well plate (Corning Cat. # 3676), in quadruplicate so rows A-D are staurosporine replicates, E-H are JAK2 Inhibitor 2 replicates, etc.

Setup Guide on the BioTek Synergy™ 2 Multi-Mode Microplate Reader

4. Add 2.5 µl of kinase buffer alone to rows 21 and 22 (0% inhibition no compound control), 23 (0% phosphorylation control, no kinase added) and 24 (Phosphopeptide 100% phosphorylation positive control).
5. Add 5 µL of the 2X Peptide/Kinase Mixture (800 nM Fluorescein-Poly GT peptide, PV3610, 222 ng/ml JAK2 JH1/JH2 or 270 ng/ml JAK2 JH1/JH2 V617F, determined experimentally as outlined above) to Columns 1-22. DO NOT ADD TO COLUMN 23 OR 24. Add 5 µL of 800 nM substrate alone without kinase to Column 23, rows A-L (0% phosphorylation control) and 5 µl of 800 nM phosphopeptide control substrate to Column 24, rows A-L (100% phosphorylation control). Add 5µl kinase buffer alone to the remaining 8 wells (Columns 23 and 24, rows M-P) as a buffer-only reference.
6. Add 2.5 µL of 4X ATP Solution (20 µM) per well to all Columns to start reaction.
7. Shake assay plate on a plate shaker for 30 seconds.
8. Incubate assay plate for 60 minutes at room temperature.
9. Add 10 µl per well of 20 mM EDTA, 2 nM TB-PY20 antibody (PV3552) mix diluted in TR-FRET Dilution Buffer (PV3574) per well to stop kinase reaction. Do not add to Rows M-P, Columns 23 and 24 (buffer controls only), instead replace with 10 µl TR-FRET Dilution Buffer supplemented with EDTA only.
10. Shake plate again on a plate shaker for 30 seconds.
11. Incubate for 30 minutes at room temperature.
12. Read and analyze as directed in the protocol.

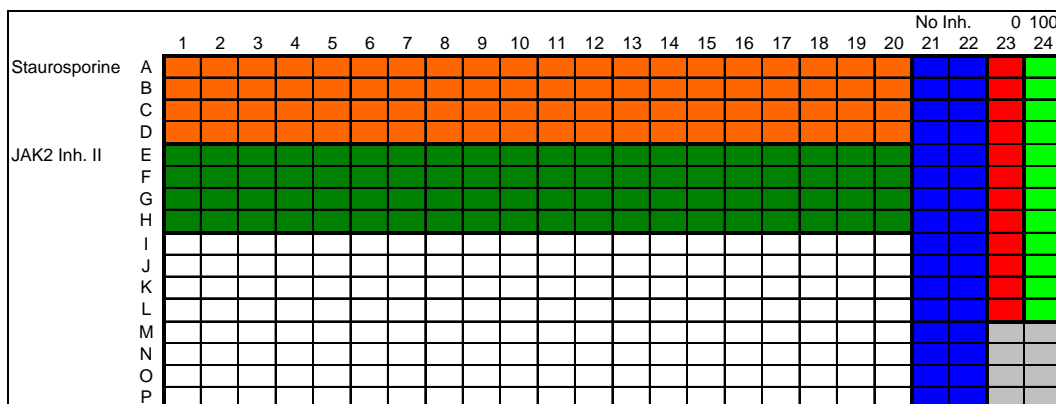


Figure 2: Assay Plate Schematic. Compound titrations shown in Columns 1-20, Columns 21 and 22 prepared without any inhibitor as kinase activity controls, Column 23 prepared with no kinase (0% phosphorylation) and Column 24 prepared using phosphopeptide control (100% phosphorylation). Note 8 wells in gray in bottom right, which were prepared without any inhibitor, substrates, or antibody as buffer controls.

D. Results

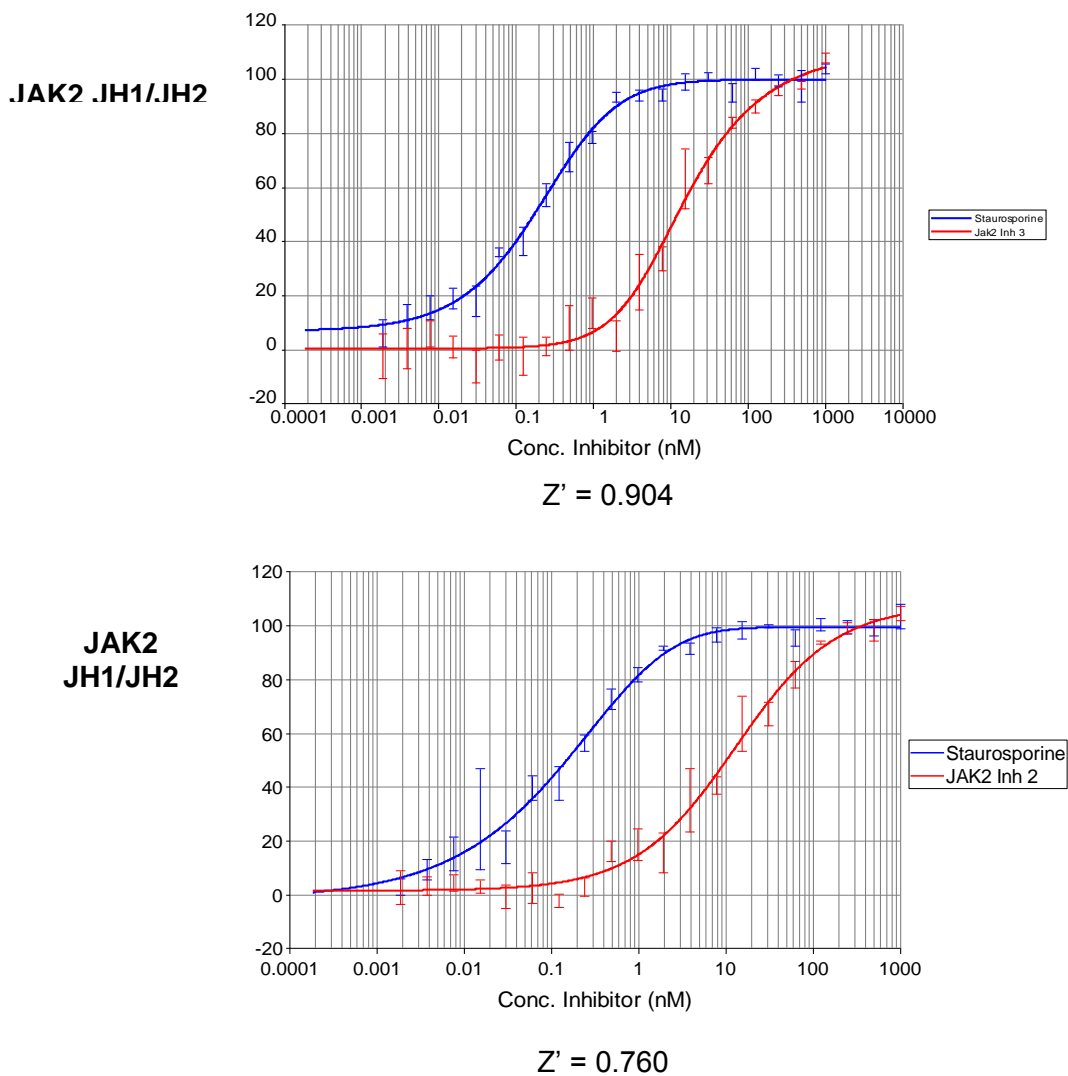


Figure 3: LanthaScreen® Assay. LanthaScreen® assay performed on the BioTek Instruments Synergy™ 2.