

## Z'-LYTE<sup>®</sup> Assay Setup Guide on the Tecan Safire<sup>2™</sup> Microplate Reader

NOTE: The Tecan Safire<sup>2™</sup> Microplate Reader was tested for compatibility with Invitrogen's Z'-LYTE<sup>®</sup> Assay using the Z'-LYTE<sup>®</sup> Tyr6 kit (PV4122) against JAK2 JH1/JH2 and JAK2 JH1/JH2 V617F kinases. The following document is intended to demonstrate setup of this instrument. For more detailed information and technical support of Invitrogen assays please call 1-800-955-6288, select option "3", then extension 40266. For more detailed information and technical support of Tecan instruments or software, please contact Tecan at 1-888-798-0538 or [info@tecan.com](mailto:info@tecan.com).

### A. Recommended Optics

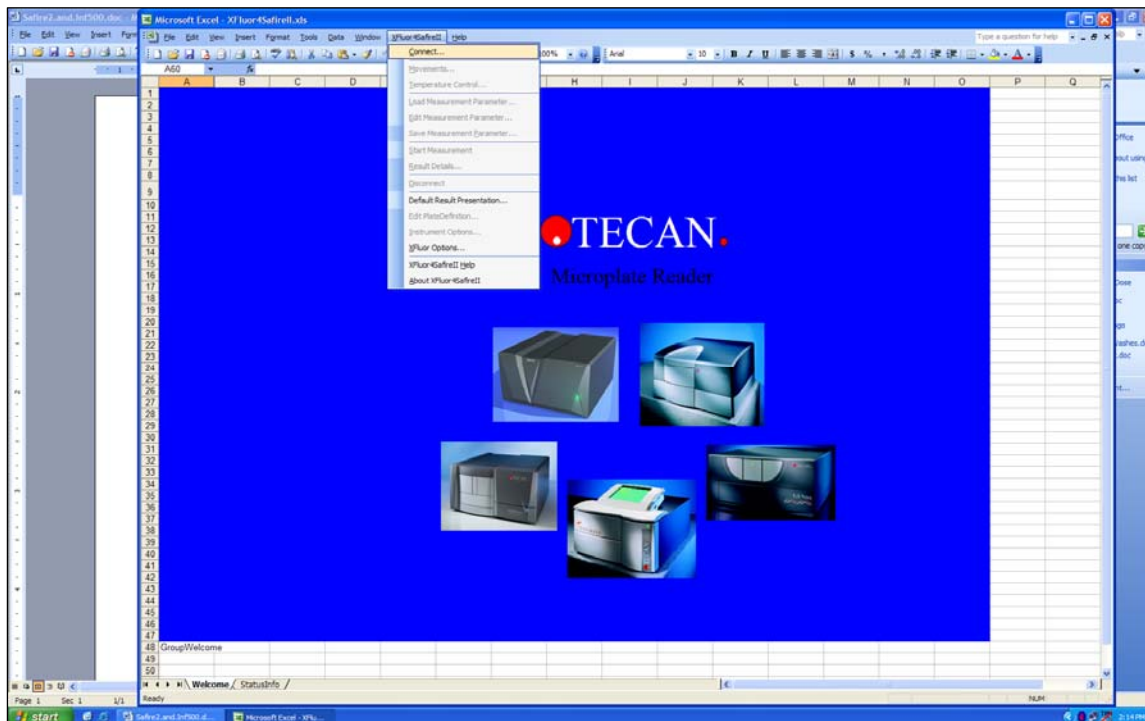
	wavelength (nm)	diameter (mm)
Excitation	405/12	monochromator
Emission 1	445/12	monochromator
Emission 2	520/12	monochromator

### B. Instrument Setup

1. Make certain plate reader is turned on, and open up XFluor Data Manager software on computer.

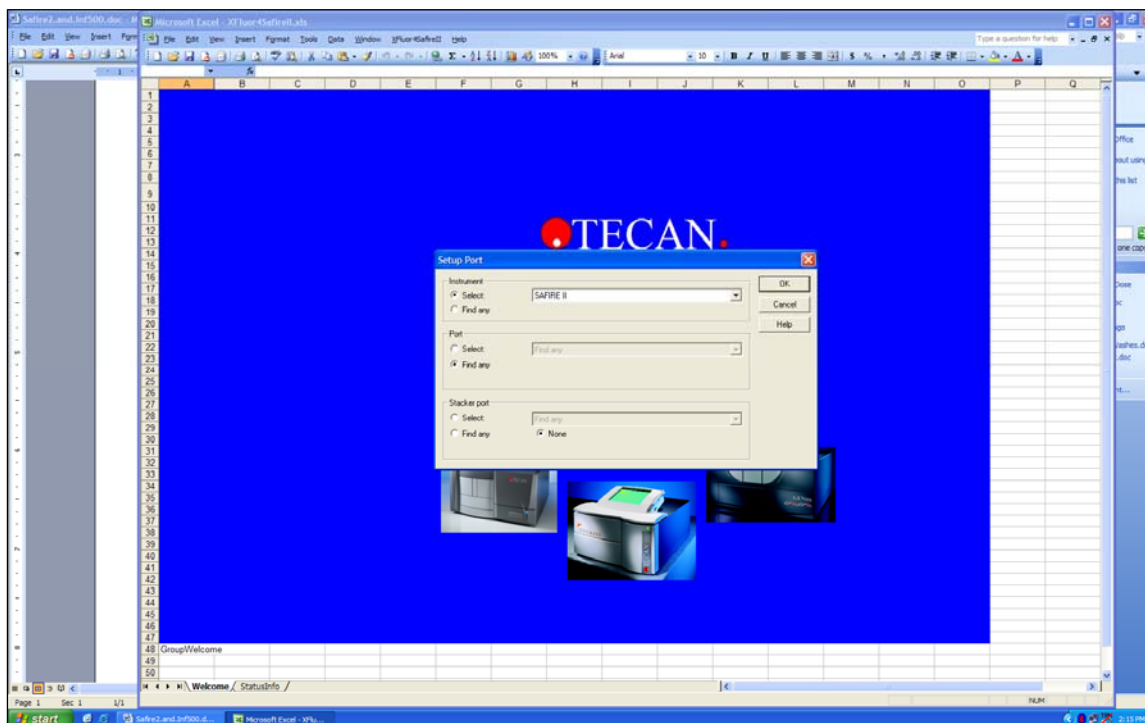
## Setup Guide on the Tecan Safire<sup>2</sup>™ Microplate Reader

2. When XFluor opens, it will default to a generic starting page. From the menu bar on the top, go to "XFluor4SafireII" and select "Connect" from the drop-down menu.



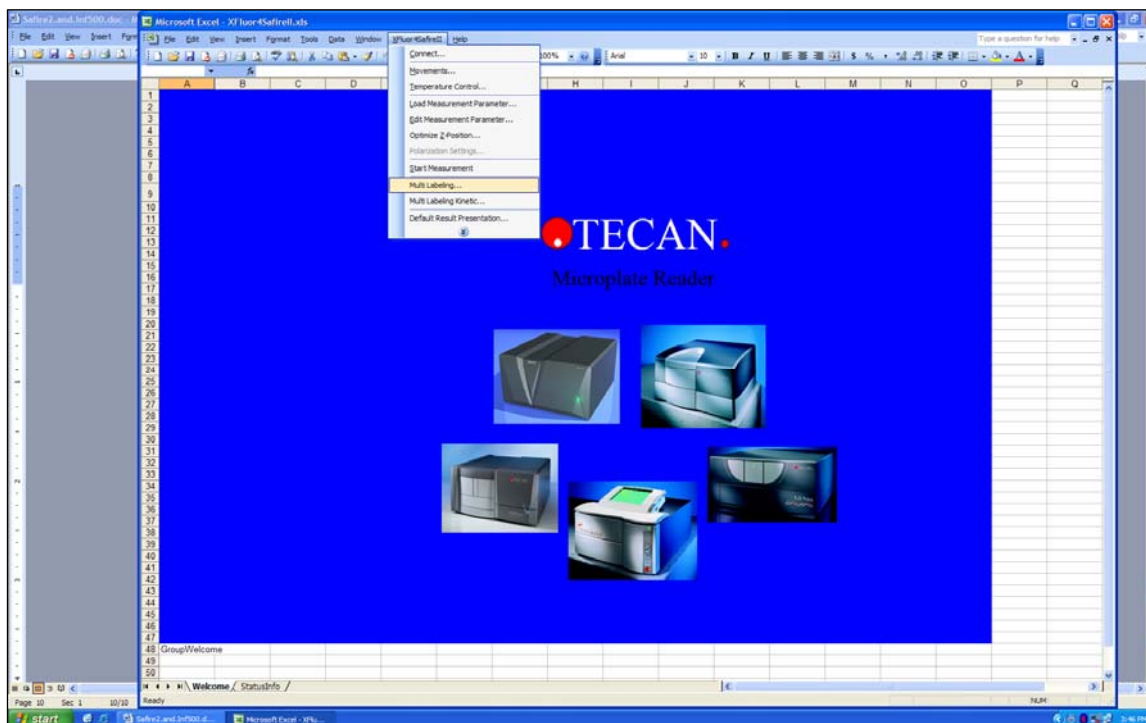
## Setup Guide on the Tecan Safire<sup>2</sup>™ Microplate Reader

3. Select your instrument (Safire II) and click OK.



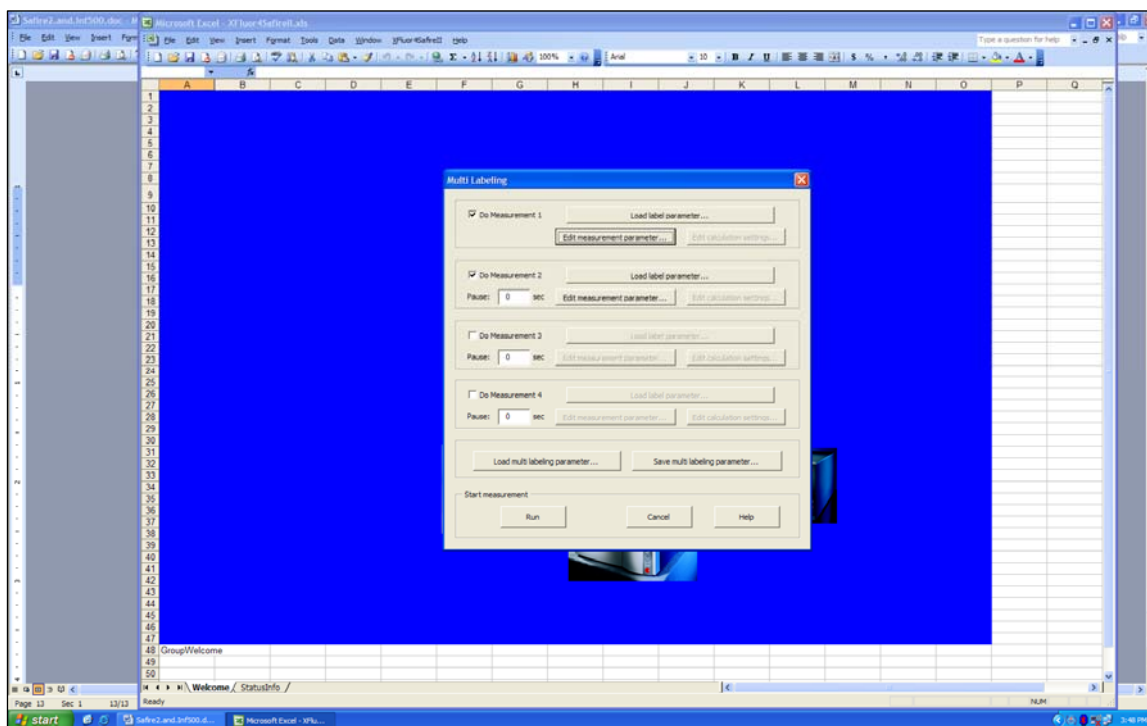
## Setup Guide on the Tecan Safire<sup>2</sup>™ Microplate Reader

4. From the XFluor drop-down menu, select "Multi Labeling".



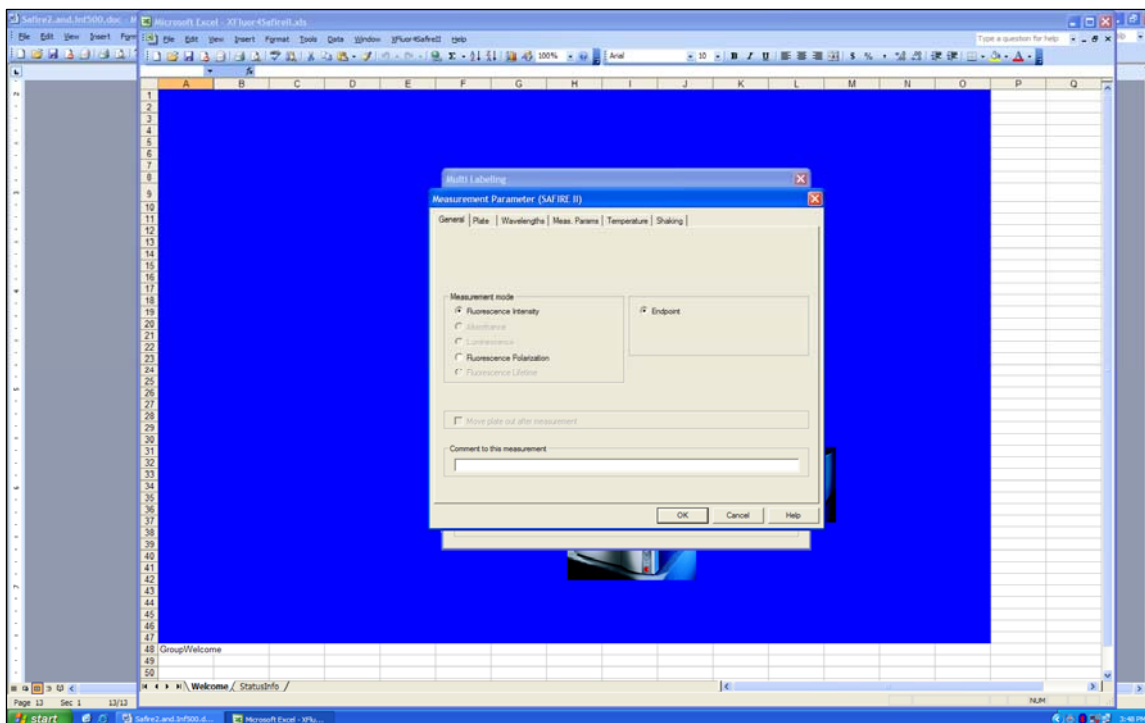
**Setup Guide on the Tecan Safire<sup>2</sup>™ Microplate Reader**

5. A new window will open. Z'-LYTE® uses a donor and an acceptor, so check the boxes for "Do Measurement 1" and "Do Measurement 2". Next, click the "Edit Measurement Parameter" tab in the Measurement 1 box to set up the donor fluor measurement parameters.



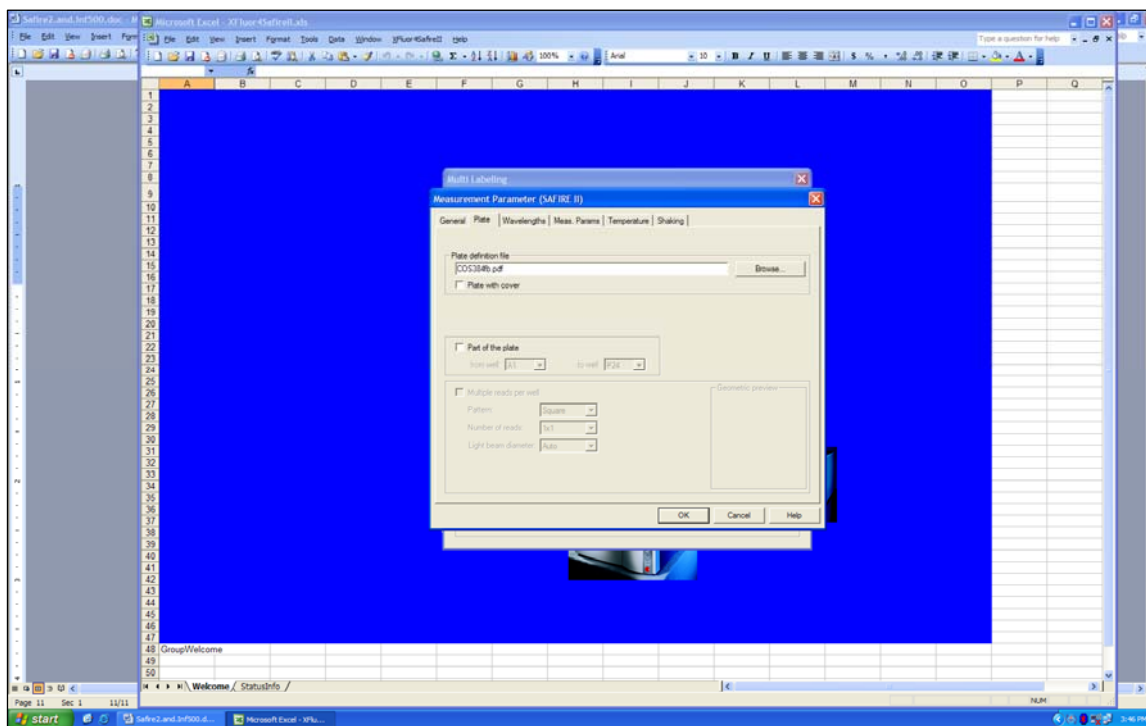
## Setup Guide on the Tecan Safire<sup>2</sup>™ Microplate Reader

6. A new window will again open; under the General tab make sure "Fluorescence Intensity" and "Endpoint" are selected.



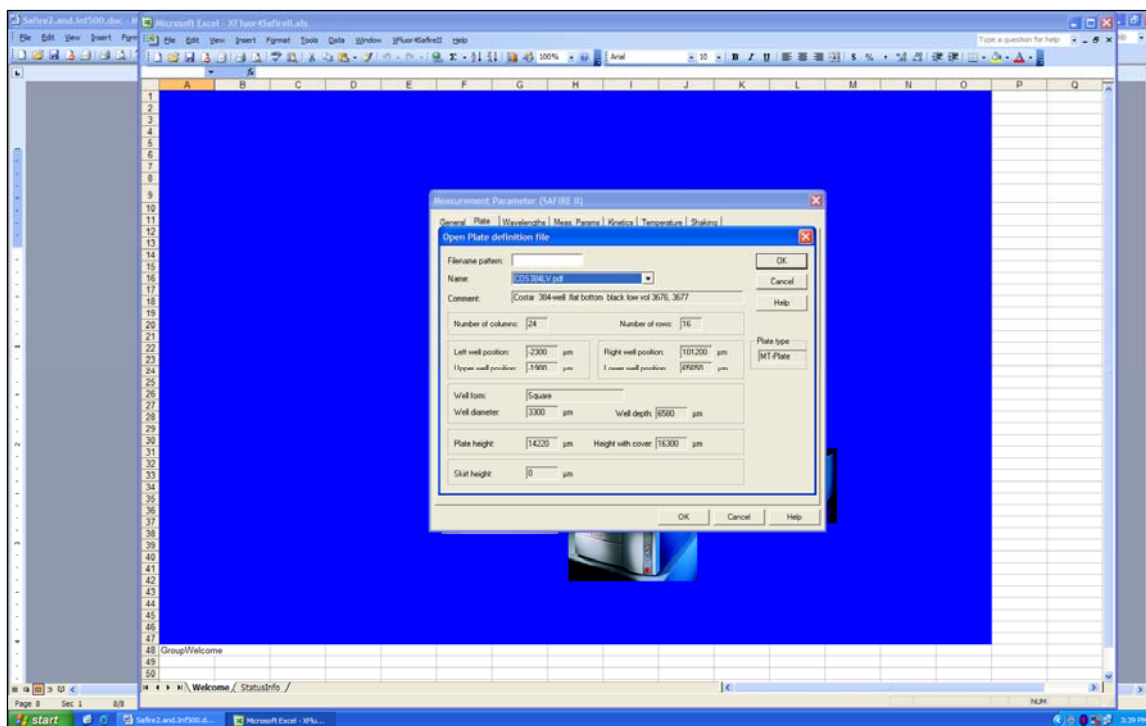
## Setup Guide on the Tecan Safire<sup>2</sup>™ Microplate Reader

7. Select the Plate tab. Make sure the whole plate is set to be read, and click the "Browse" tab to select your plate.



## Setup Guide on the Tecan Safire<sup>2</sup>™ Microplate Reader

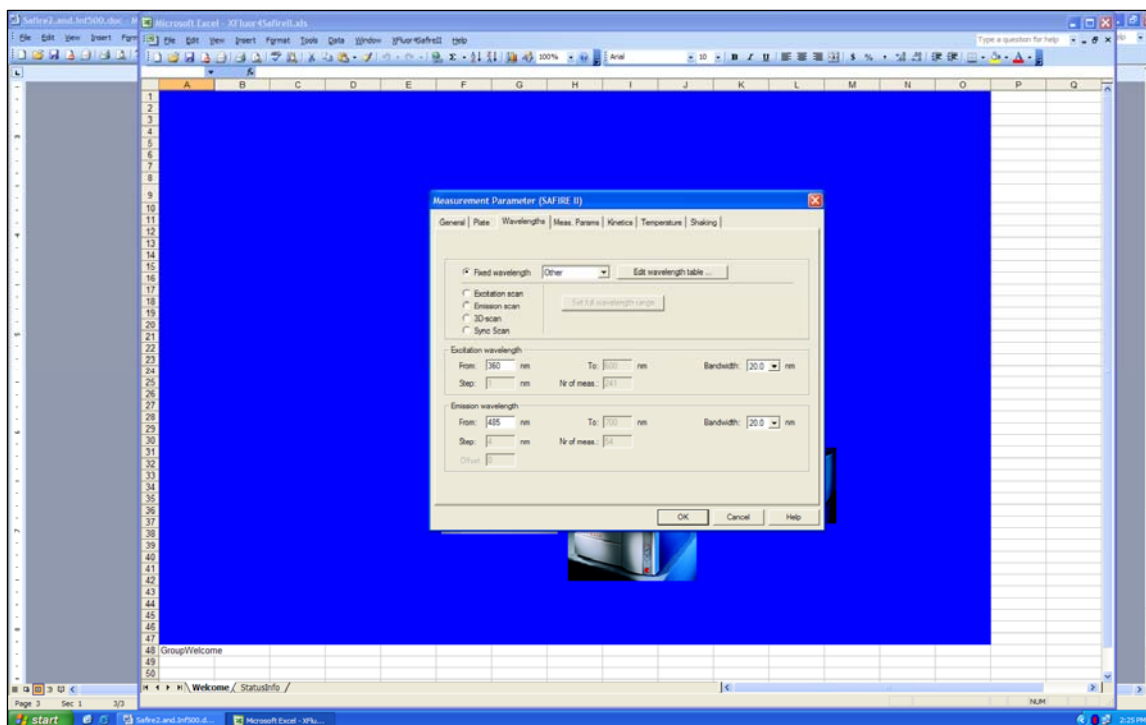
8. A new window will appear; from the drop-down menu, select your plate. When finished, select OK to close this window.





**Setup Guide on the Tecan Safire<sup>2</sup>™ Microplate Reader**

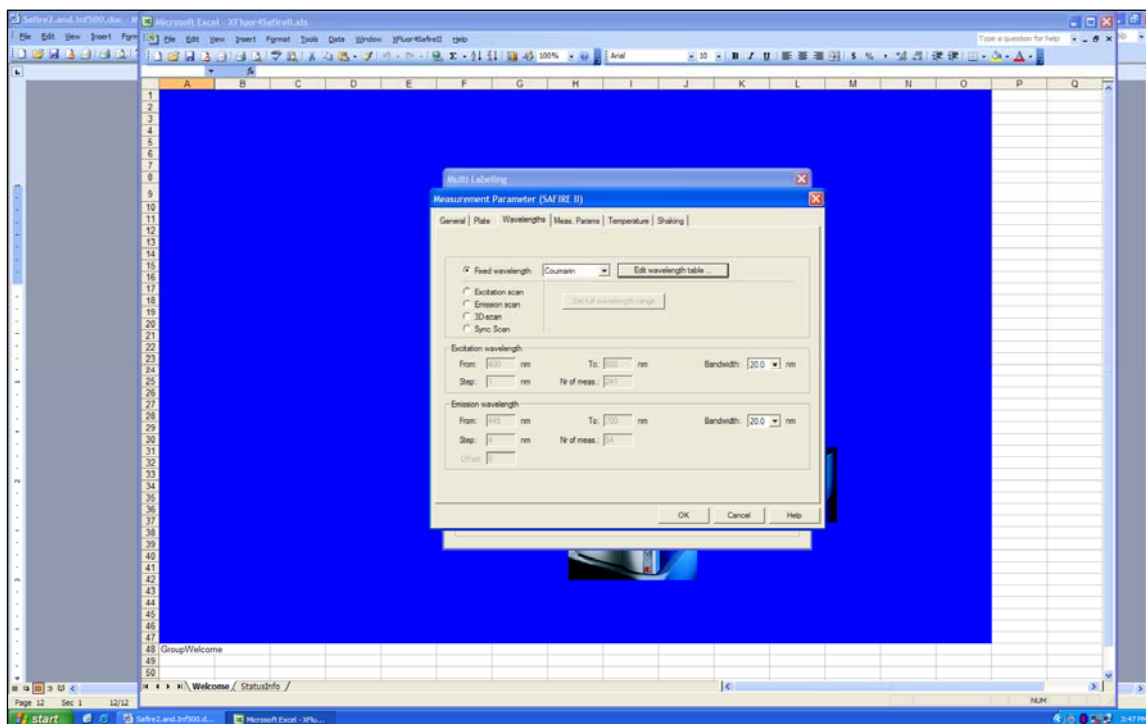
9. Select the Wavelengths tab. Check the "Fixed Wavelength" button, and then select the "Edit wavelength table" tab to enter your excitation and donor emission settings.





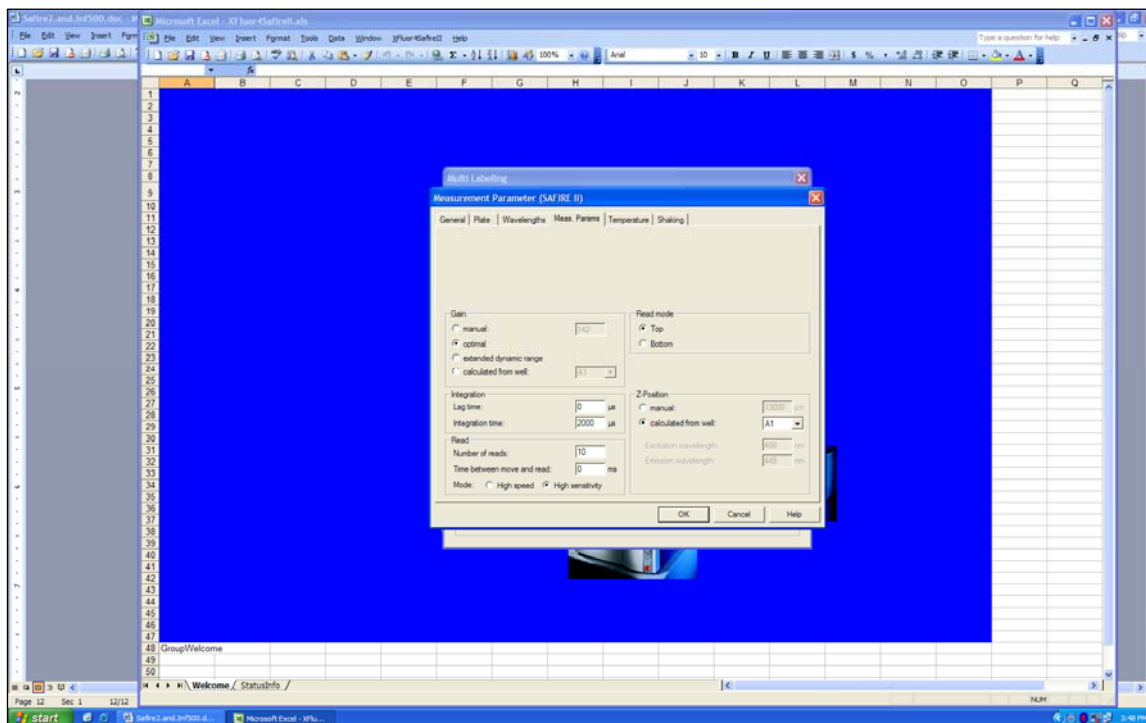
## Setup Guide on the Tecan Safire<sup>2</sup>™ Microplate Reader

11. You will return to the Measurement Parameter window. Select the appropriate bandwidth settings and when finished select the Meas. Params. tab.



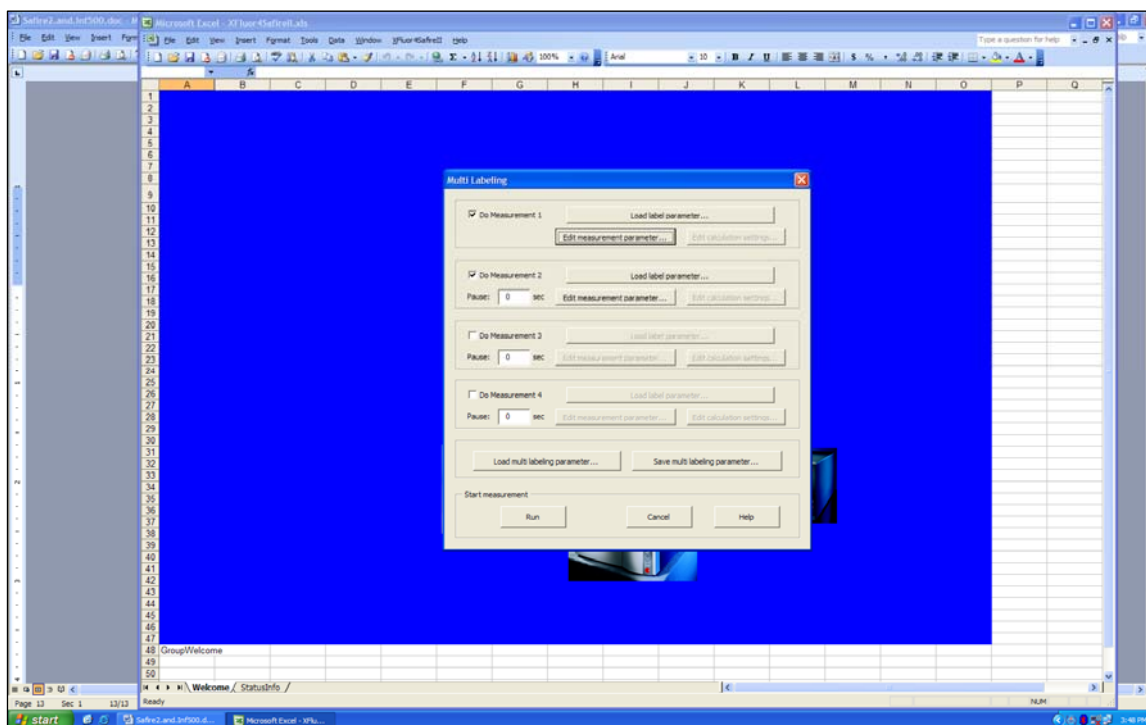
**Setup Guide on the Tecan Safire<sup>2</sup>™ Microplate Reader**

12. Make sure the Read mode is set to Top, and in this case Gain was set to Optimal and Z-Position was set to well A1. When finished with all parameters, click OK at the bottom right.



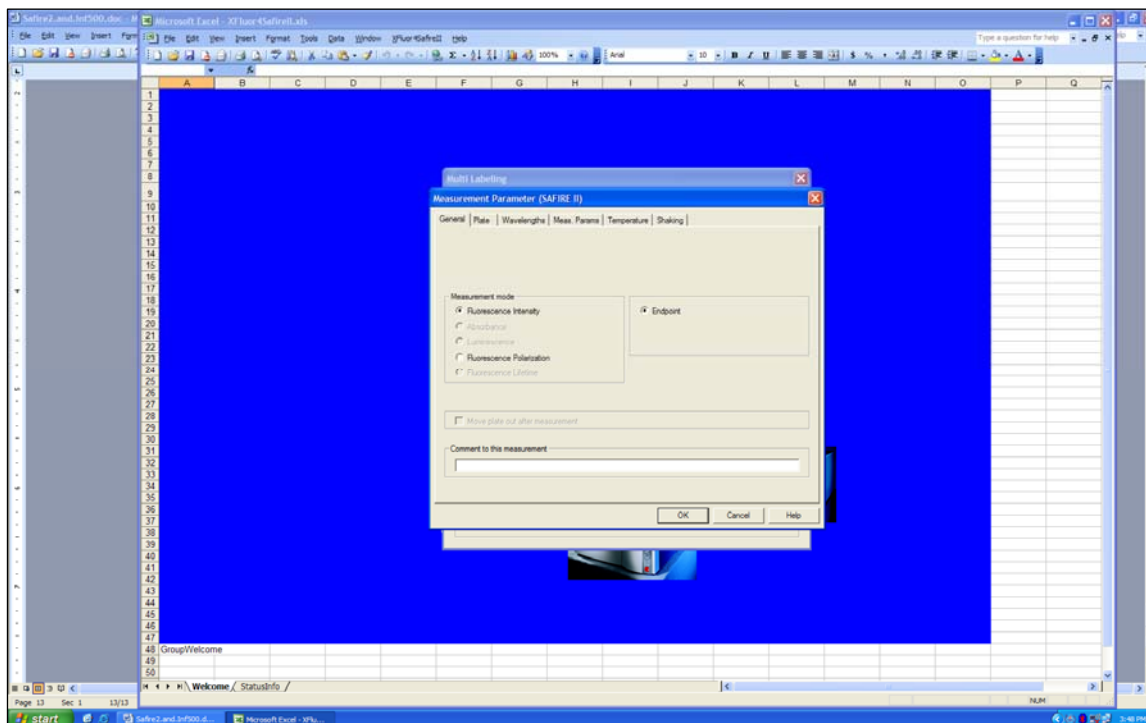
**Setup Guide on the Tecan Safire<sup>2</sup>™ Microplate Reader**

13. XFluor will default back to the Multi Labeling window. Select the "Edit Measurement Parameter" tab beneath the checked Do Measurement 2 box to edit the acceptor fluor measurement parameters.



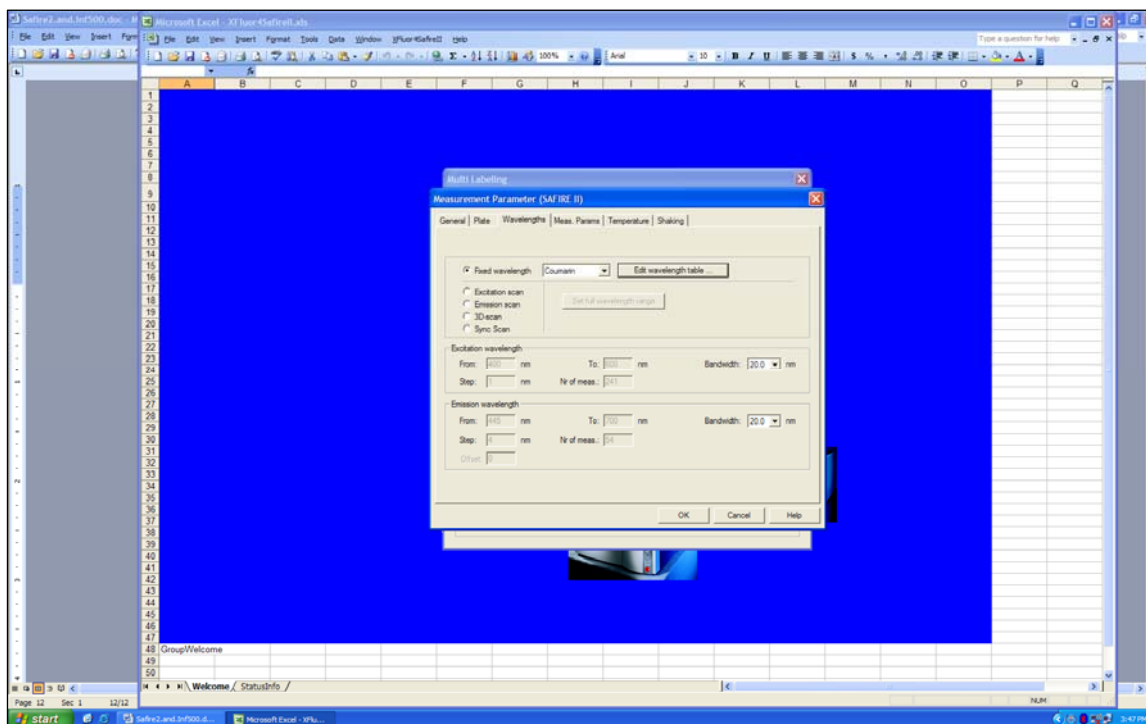
**Setup Guide on the Tecan Safire<sup>2</sup>™ Microplate Reader**

14. Under the General tab, make sure "Fluorescence Intensity" and "Wavelengths" are checked. At this point the plate window will be defaulted because of the first set of donor parameters, so skip the Plate tab and open the Wavelengths tab.



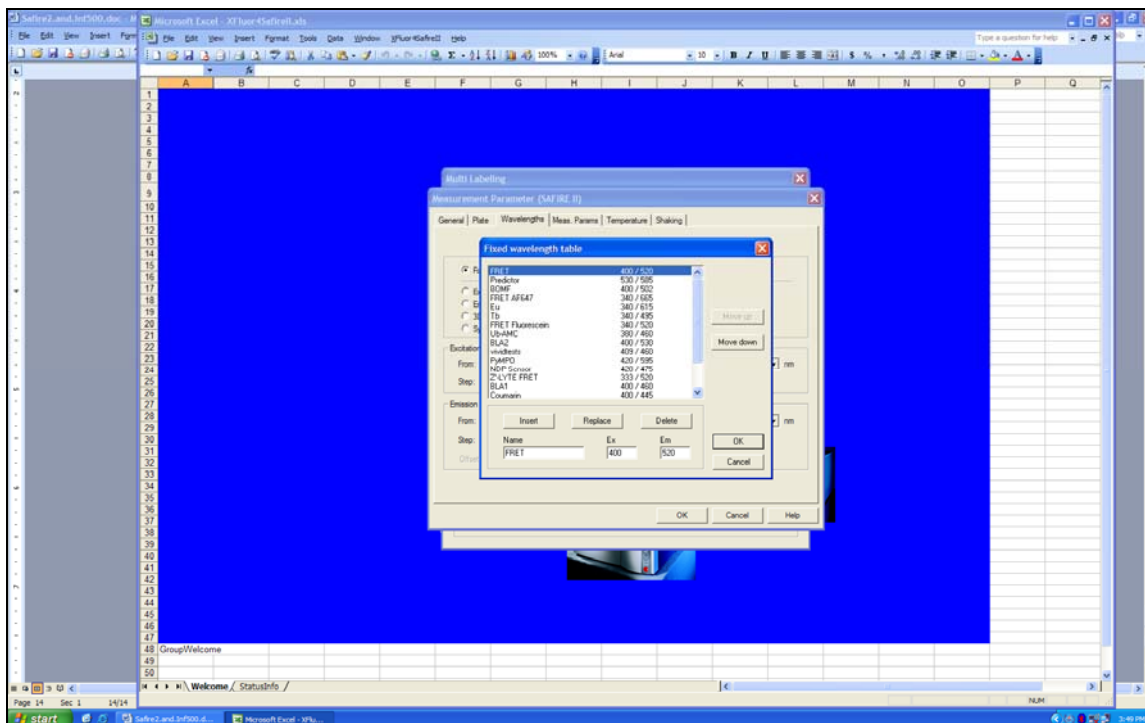
## Setup Guide on the Tecan Safire<sup>2</sup>™ Microplate Reader

15. As before, select the "Edit wavelength table" tab.



## Setup Guide on the Tecan Safire<sup>2</sup>™ Microplate Reader

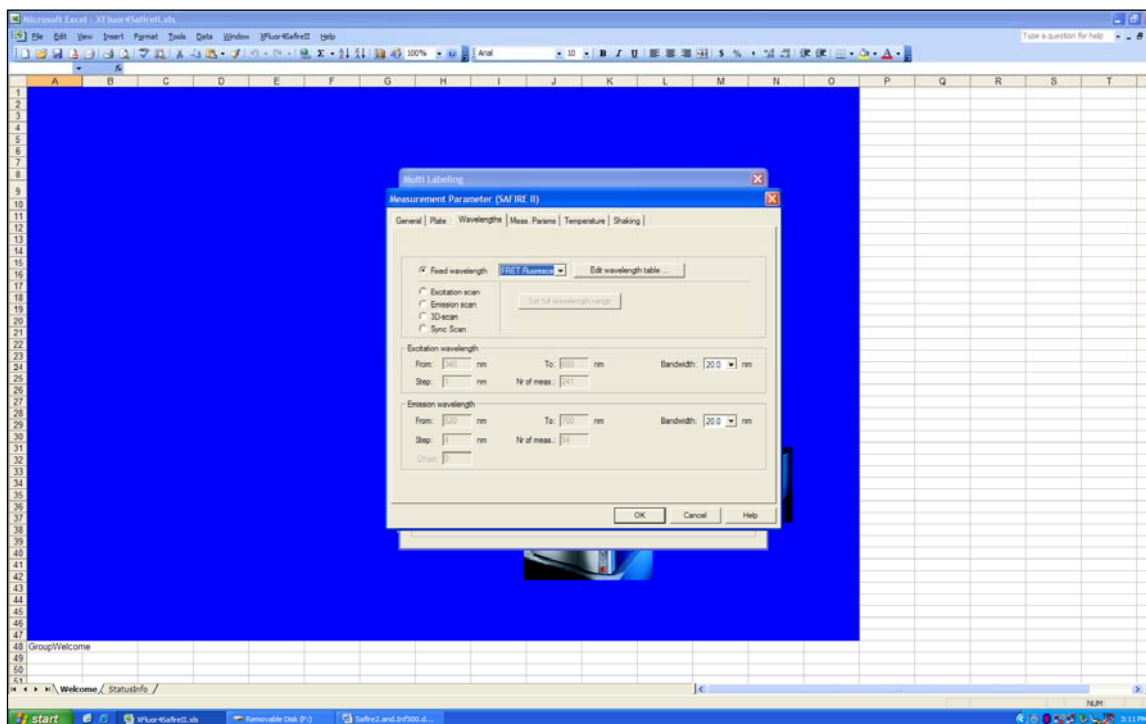
16. In the popup Fixed wavelength table enter a name and excitation and emission values for the Z'-LYTE<sup>®</sup> acceptor. When finished select "Insert" to add them to the table and then select OK.





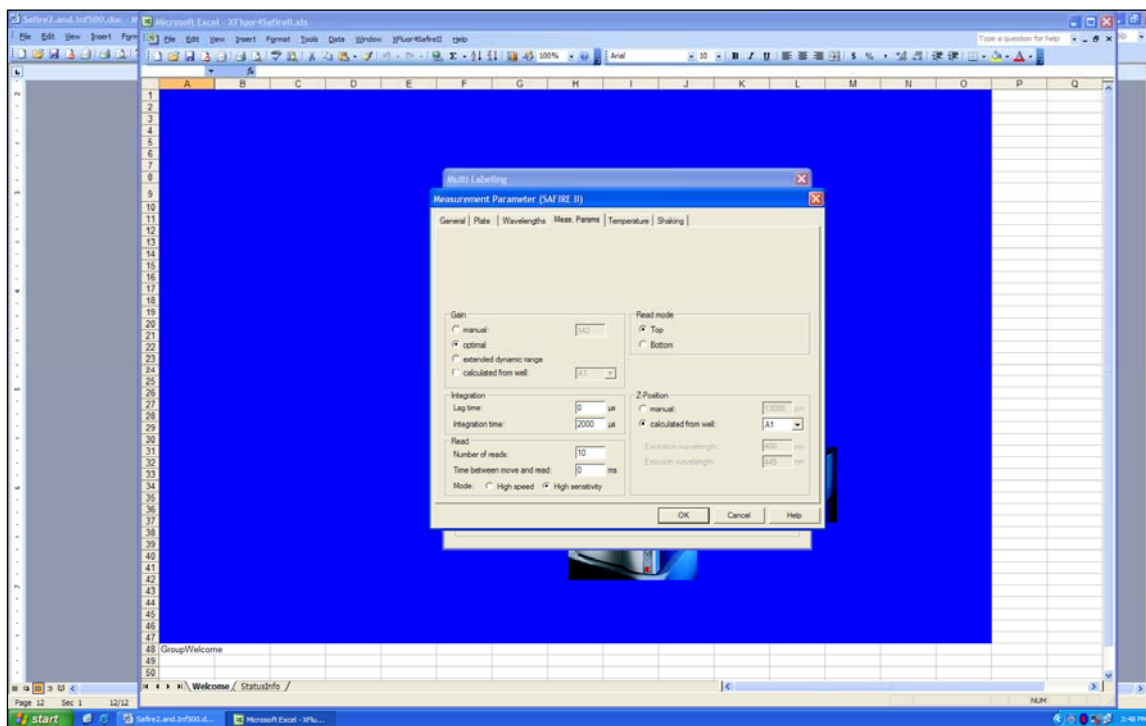
## Setup Guide on the Tecan Safire<sup>2</sup>™ Microplate Reader

17. Back in the Measurement Parameter window, select bandwidths and then select the Meas. Params tab.



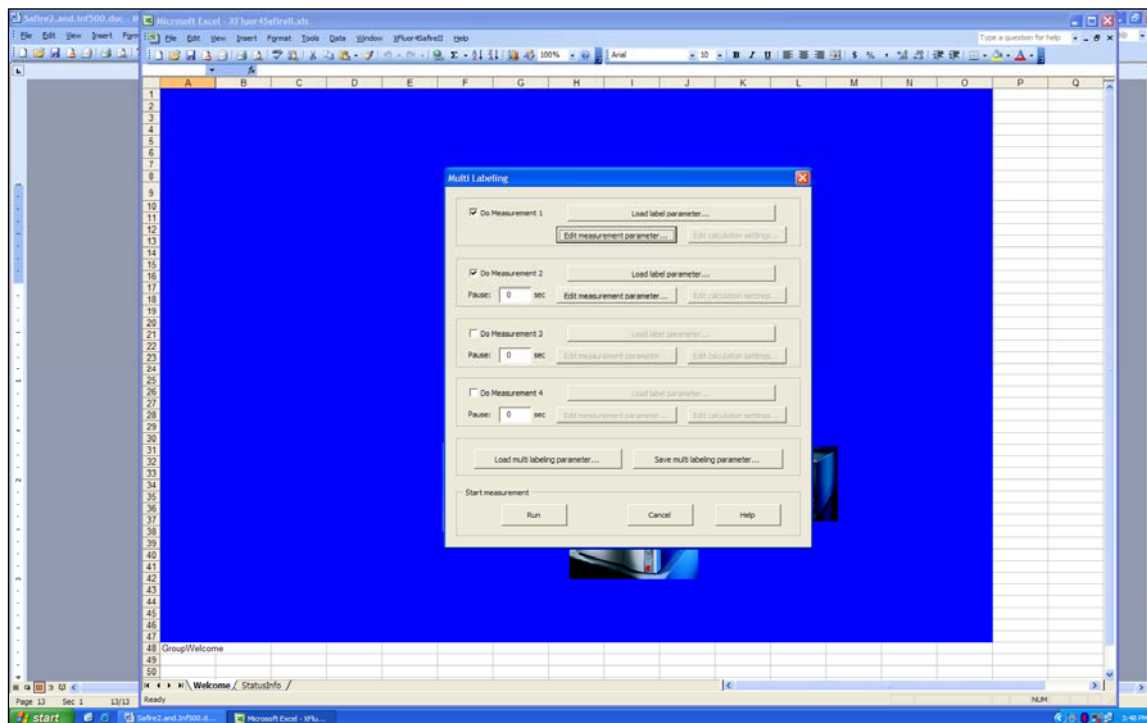
## Setup Guide on the Tecan Safire<sup>2</sup>™ Microplate Reader

18. Set the Gain and Z-Position as shown below. Make sure Read mode is set to "Top". When finished, select OK.



**Setup Guide on the Tecan Safire<sup>2</sup>™ Microplate Reader**

19. Once again the program will default back to the initial Multi Labeling window. Make sure plate is inserted (or insert now, in "Movements" under the XFluor drop-down menu) and select "Run" to read assay plate.



## C. Z'-LYTE® Kinase Assay using JAK2 JH1/JH2 and JAK2 JH1/JH2 V617F

NOTE: The following is a sample assay performed for demonstration purposes. The section below describes how the data was obtained, and is not intended for use as an assay protocol. We recommend all first-time users follow the appropriate protocols and/or validation packets provided with their specific assay kits, and include all proper controls. The instrument settings above would be sufficient for any Z'-LYTE® assay, the information below is provided as representative data. Assay was run at ATP  $K_{mapparent}$  and a kinase concentration producing approximately 30-40% of maximal phosphorylation, as discussed in Section 9 and 10 of the Z'-LYTE® protocols. ATP and kinase concentrations should be optimized for each kinase by the actual user. Specific Z'-LYTE® assay protocols and setup information from Invitrogen's own in-house SelectScreen® Custom Profiling Z'-LYTE®-based kinase assay service can be located at the following link: <http://www.invitrogen.com/content.cfm?pageid=9866>.

1. Prepare initial 100X serial dilution curves in rows A and E of a 384-well plate: Dilute Staurosporine and JAK2 Inhibitor II to a 100X initial concentration in 100% DMSO (100  $\mu$ M). Prepare a set of 1:1 serial dilutions from the initial concentration in a 384-well plate, starting with 80  $\mu$ l in Column 1 and 40  $\mu$ l DMSO in wells 2-20. Add 40  $\mu$ l from well 1 to well 2, and then mix well 2, and take 40  $\mu$ 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
Staurosporine	100X	A																							
	4X	B																							
		C																							
		D																							
JAK2 Inh. II		E																							
	4X	F																							
		G																							
		H																							
		I																							
		J																							
		K																							
		L																							
		M																							
		N																							
		O																							
		P																							

**Figure 1: Schematic of initial compound dilution.** Staurosporine and JAK2 Inhibitor II were titrated from a 100  $\mu$ 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 (50 mM HEPES pH 7.5, 0.01% BRIJ-35, 10 mM MgCl<sub>2</sub>, 1 mM EGTA) in the row below by adding 2  $\mu$ l of diluted inhibitor from the well above to

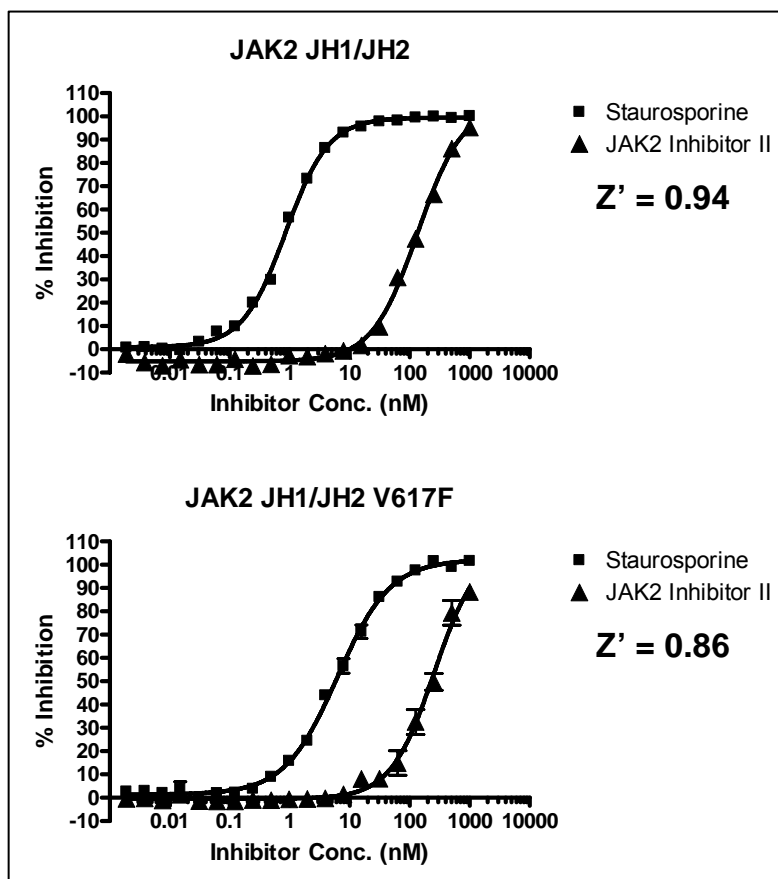
- 48 µl of kinase buffer. This will produce a final serial dilution starting at 4 µM, which will then produce a final assay concentration starting at 1 µM.
3. Begin to prepare an assay plate: Add 2.5 µL of the compound dilutions per well into a low volume NBS, 384-well plate (Corning Cat. # 3676), in quadruplicate so rows A-D are staurosporine replicates, E-H are JAK2 Inhibitor 2 replicates, etc.
  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 (2 µM Tyr 06 peptide, 2600 ng/ml JAK2 JH1/JH2 or 1300 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 2µM substrate alone without kinase to Column 23, rows A-L (0% phosphorylation control) and 5 µl of 2 µM phpsphopeptide control substrate to Column 24, rows A-L (100% phosphorylation control). Add 5 µl 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 (200 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 5 µL of the Development Reagent Solution to each well. Use the lot-specific dilutions indicated on your CoA as dilution may vary based upon Z'-LYTE® peptide and Development Reagent A lot.
  10. Shake plate again on a plate shaker for 30 seconds.
  11. Incubate for 60 minutes at room temperature.
  12. Read and analyze as directed in the protocol.

### Setup Guide on the Tecan Safire<sup>2</sup>™ Microplate Reader

		1	2	3	4	5	6	7	8	9	10	11	12	13	14	15	16	17	18	19	20	No Inh.		0	100
Staurosporine	A																								
	B																								
	C																								
	D																								
JAK2 Inh. II	E																								
	F																								
	G																								
	H																								
	I																								
	J																								
	K																								
	L																								
	M																								
	N																								
	O																								
	P																								

**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 with out any inhibitor or substrates, as buffer controls.

## D. Results



**Figure 1: Z'-LYTE® Kinase Assay.** Z'-LYTE® assay performed using the Tecan Safire<sup>2</sup>™.