

## GeneBLazer® Assay Setup Guide on the BMG LABTECH OPTIMA Microplate Reader

NOTE: The BMG LABTECH OPTIMA Microplate Reader was tested for compatibility with Invitrogen's GeneBLazer® Assay in bottom-read mode using two CellSensor® cell lines; the *irf1-bla* HEL and *irf1-bla* TF-1 cell lines (K1647 and K1657, respectively). The following document is intended to demonstrate setup of this instrument and provide representative data. 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 BMG LABTECH instruments or software, please contact BMG LABTECH at 1-877-264-5227 or [www.bmglabtech.com](http://www.bmglabtech.com).

### A. Recommended Optics

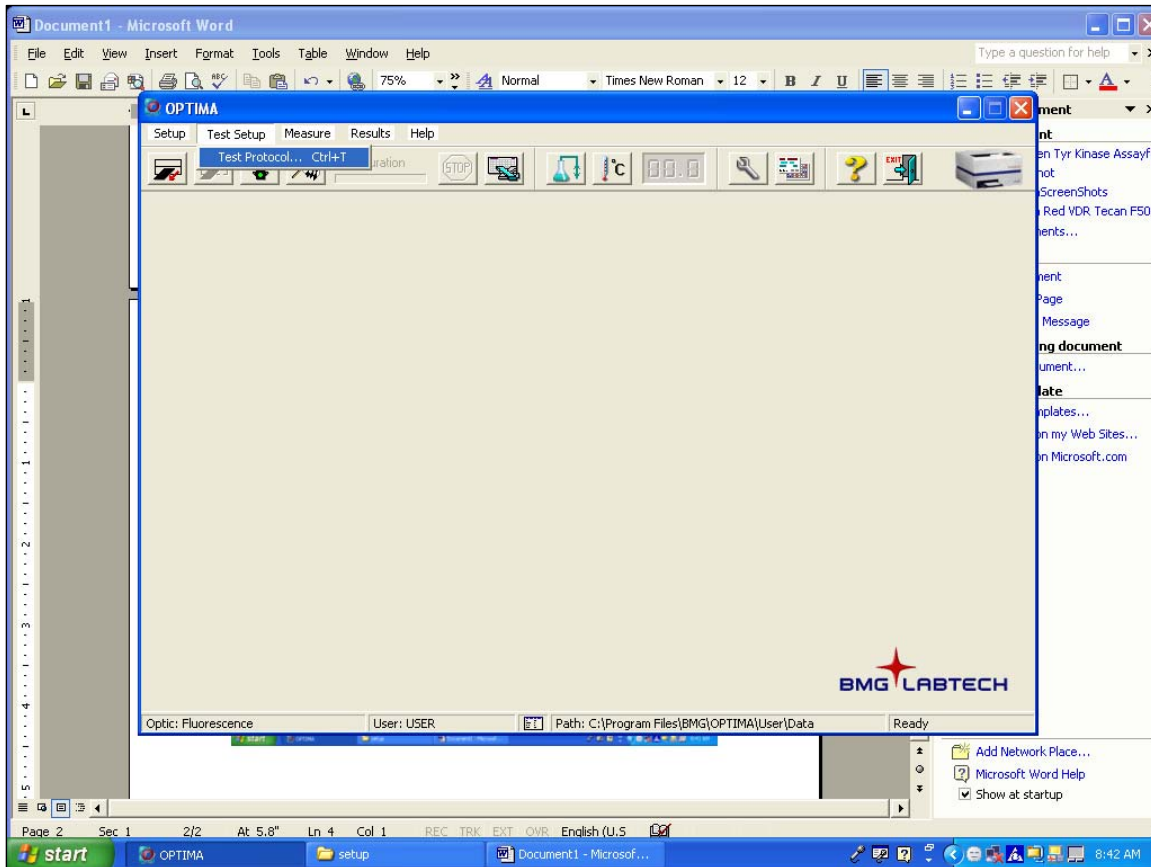
	wavelength (nm)	BMG LABTECH Filters
Excitation	400 (or similar)	*contact BMG LABTECH
Emission 1	440 (or similar)	*contact BMG LABTECH
Emission 2	520 (or similar)	*contact BMG LABTECH

### B. Instrument Setup

1. Make certain plate reader is turned on, and open up OPTIMA Control software on computer. Insert plate into plate reader.

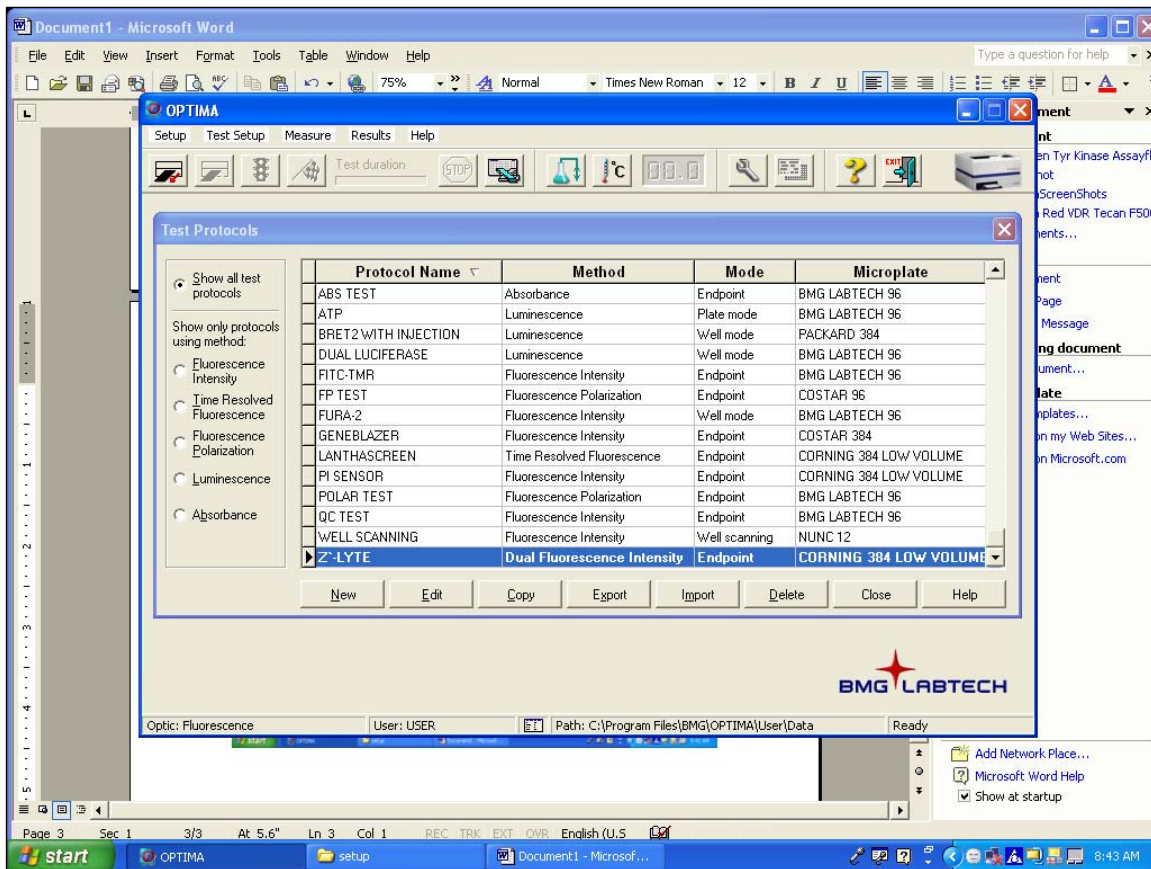
**Setup Guide on the BMGLABTECH OPTIMA Microplate Reader**

- When OPTIMA Control software opens, if you do not have a pre-existing protocol for GeneBLAzer®, select "Test Protocol" from the Test Setup menu bar at the top of the window.



**Setup Guide on the BMGLABTECH OPTIMA Microplate Reader**

- At this point, a new screen will open (below). Click on the “Show all test protocols” button on the left side of the screen, then select “New” from the tabs at the bottom.

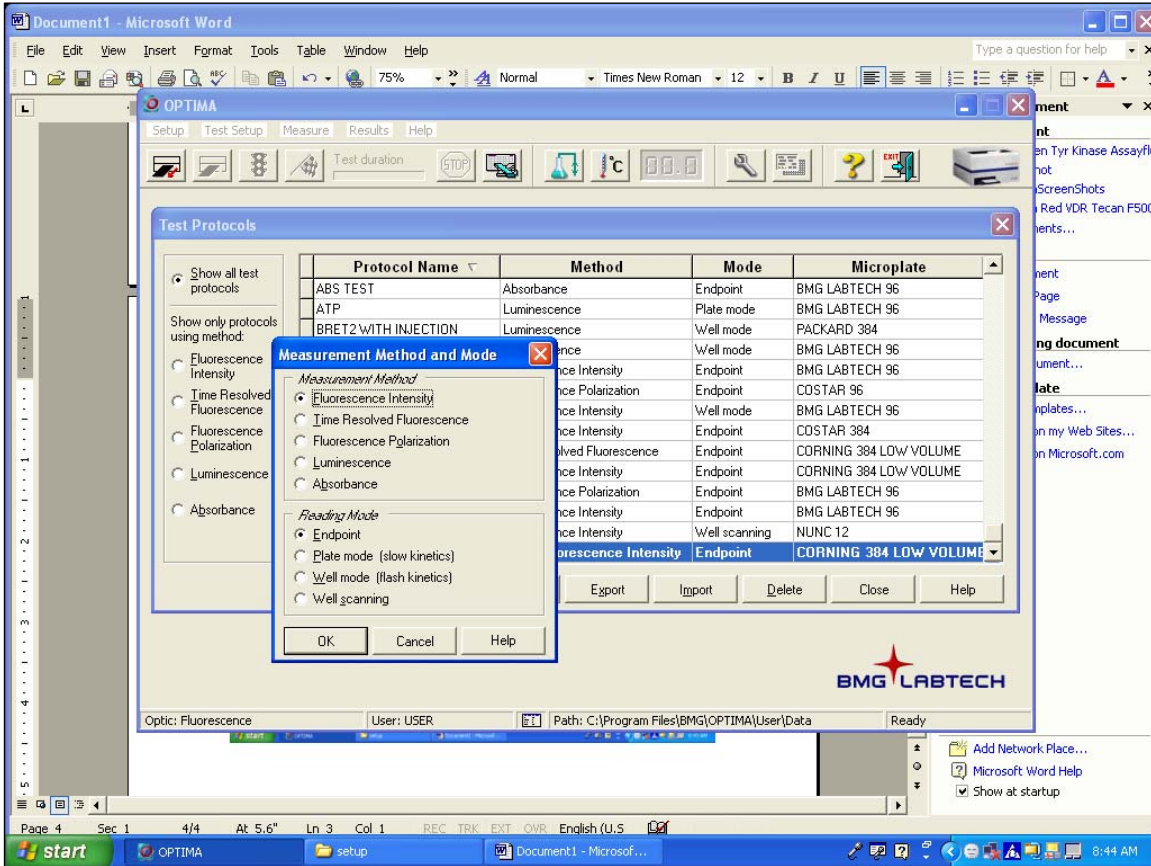


Have a question? Contact our Technical Support Team

NA: 800-955-6288 or INTL: 760-603-7200 Select option 3, ext. 40266 Email: drugdiscoverytech@invitrogen.com

**Setup Guide on the BMGLABTECH OPTIMA Microplate Reader**

4. A new window will pop up. Select “Fluorescence Intensity” and “Endpoint” and then select “OK”.

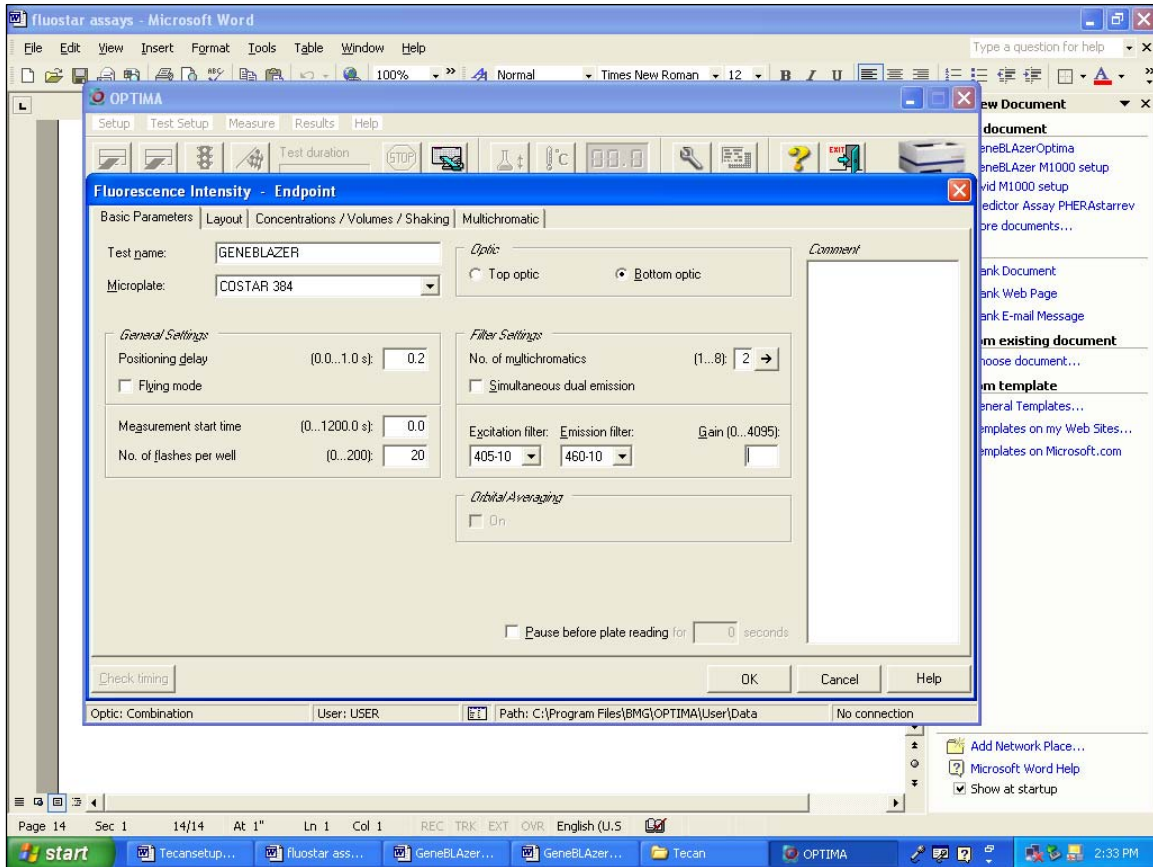


Have a question? Contact our Technical Support Team

NA: 800-955-6288 or INTL: 760-603-7200 Select option 3, ext. 40266 Email: drugdiscoverytech@invitrogen.com

**Setup Guide on the BMGLABTECH OPTIMA Microplate Reader**

5. A new Protocol window will open automatically. Enter a test name, select plate type (there is no specification for clear-bottom plates for the equivalent 384-well plates, this does not matter), and select the **Optic (bottom)** from the drop-down list. Next, set the number of flashes per well and set up the excitation and first emission filter from the drop-down tabs. Finally, under "Filter Settings", enter "2" for the No. of multichromatics. This will allow you to click on the small arrow beside the box.



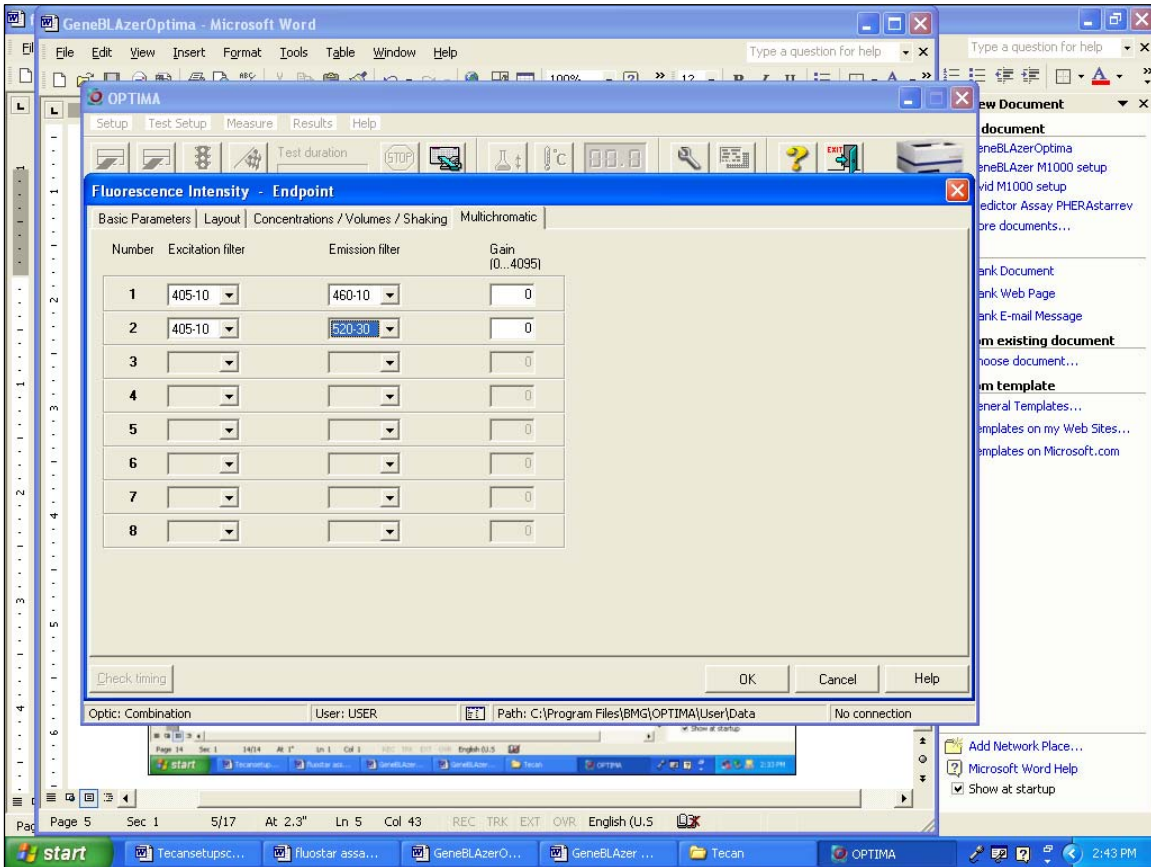
Have a question? Contact our Technical Support Team

NA: 800-955-6288 or INTL: 760-603-7200 Select option 3, ext. 40266 Email: drugdiscoverytech@invitrogen.com



**Setup Guide on the BMGLABTECH OPTIMA Microplate Reader**

- Clicking the arrow mentioned in 5 will open another window--select your filters here as shown below, then click "OK" when finished.

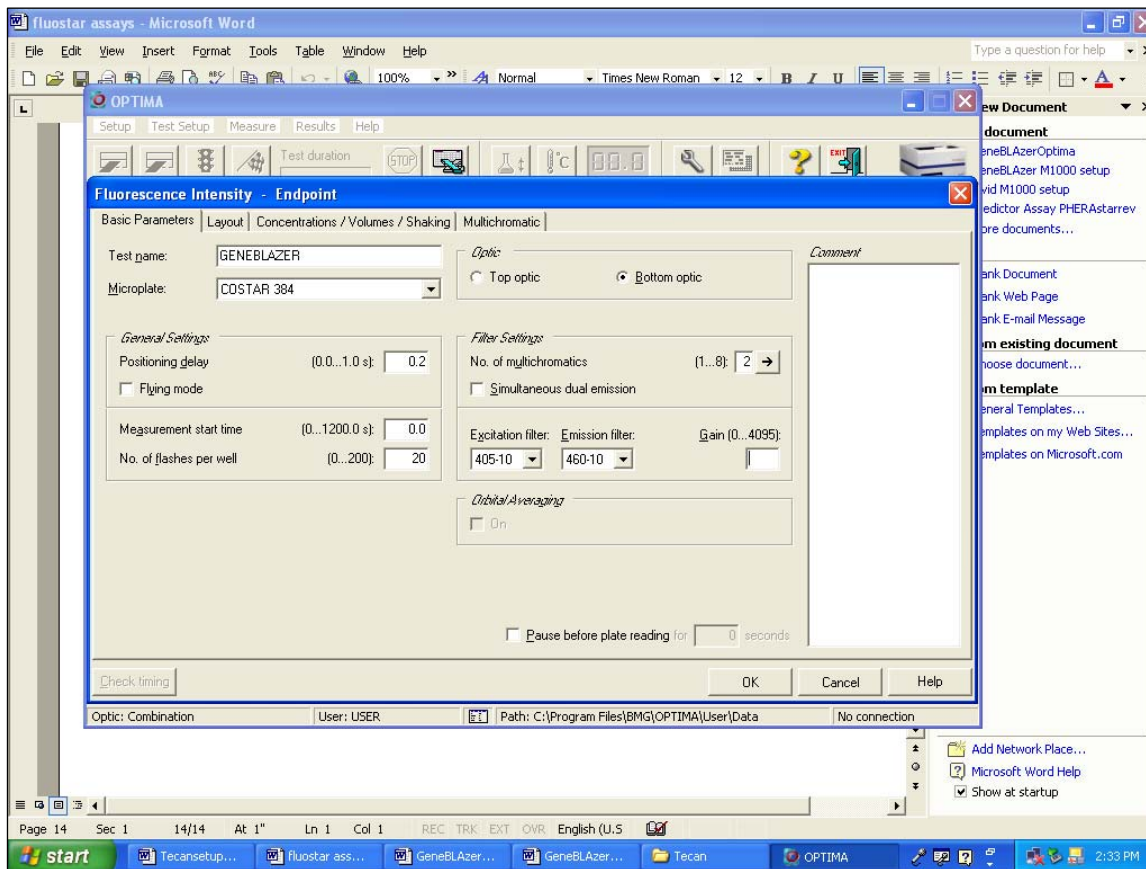


Have a question? Contact our Technical Support Team

NA: 800-955-6288 or INTL: 760-603-7200 Select option 3, ext. 40266 Email: drugdiscoverytech@invitrogen.com

**Setup Guide on the BMGLABTECH OPTIMA Microplate Reader**

7. You will return to the initial settings window. From the tabs along the top of this window, select "Layout".

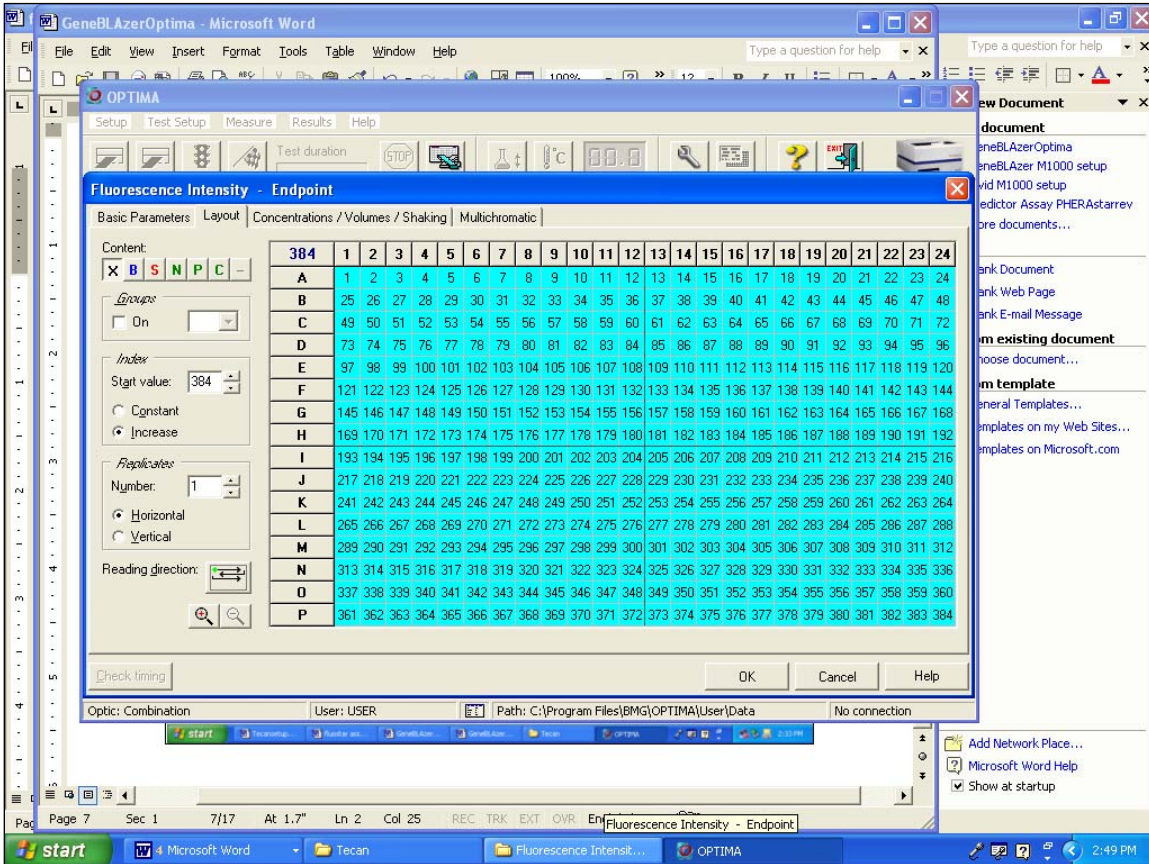


Have a question? Contact our Technical Support Team

NA: 800-955-6288 or INTL: 760-603-7200 Select option 3, ext. 40266 Email: drugdiscoverytech@invitrogen.com

**Setup Guide on the BMGLABTECH OPTIMA Microplate Reader**

- A new window will appear. Select the wells you wish to read by highlighting them. When finished, select "OK"



Have a question? Contact our Technical Support Team

NA: 800-955-6288 or INTL: 760-603-7200 Select option 3, ext. 40266 Email: drugdiscoverytech@invitrogen.com





**Figure 1: Schematic of initial compound dilution.** Staurosporine and JAK2 Inhibitor II were titrated from a 100  $\mu\text{M}$  starting concentration in the initial dilution series by preparing a 1:1 dilution curve in DMSO. A secondary dilution to 10X was then prepared in the rows below the initial dilution curve (lighter gray) using each line's specified Assay Buffer.

4. Add 36  $\mu\text{l}$  per well of cell suspension to all wells except Column 23 and 24, Rows M through P. For Column 23 and Column 24, Rows M through P, add 40  $\mu\text{l}$  Assay media alone. It is important to use cell-free wells in GeneBLAzer to background subtract donor and acceptor values for optimal GeneBLAzer assay performance.
5. Now add 4  $\mu\text{l}$  inhibitor from secondary dilutions above to wells in Columns 1-20 in quadruplicate (i.e. 4  $\mu\text{l}$  of the 10  $\mu\text{M}$  staurosporine into Column 1 Rows A through D and 4  $\mu\text{l}$  of 5  $\mu\text{M}$  staurosporine from the next well to Column 2 Rows A through D and so on to set up an actual titration from 1  $\mu\text{M}$  final starting concentration of staurosporine). Add 4  $\mu\text{l}$  Assay Buffer to Columns 21 and 22, as well as rows A-L of Columns 23 and 24, so each well has a final volume of 40  $\mu\text{l}$ .
6. Incubate HEL assay plate for 15 hours.
7. The following morning, prepare an assay plate as above for the TF-1 line, except that cells are washed once more and resuspended at  $1.5625 \times 10^6$  cells/ml in Assay Media. Add 32  $\mu\text{l}$  cells per well to all wells of the plate except Column 23 and 24, Rows M through P. For Column 23 and Column 24, Rows M through P, add 36  $\mu\text{l}$  Assay media alone, then add 4  $\mu\text{l}$  Assay Buffer containing 0.8 ng/ml GM-CSF (experimentally determined, see protocol) to all wells except Column 23 (unstimulated control) to stimulate. Add 4  $\mu\text{l}$  Assay Buffer to Column 23 to make up for the GM-CSF added to other wells. Now add 4  $\mu\text{l}$  inhibitor from secondary dilutions above to wells in Columns 1-20 in quadruplicate (i.e. 4  $\mu\text{l}$  of the 10  $\mu\text{M}$  staurosporine into Column 1 Rows A through D and 4  $\mu\text{l}$  of 5  $\mu\text{M}$  staurosporine

---

**Setup Guide on the BMGLABTECH OPTIMA Microplate Reader**

---

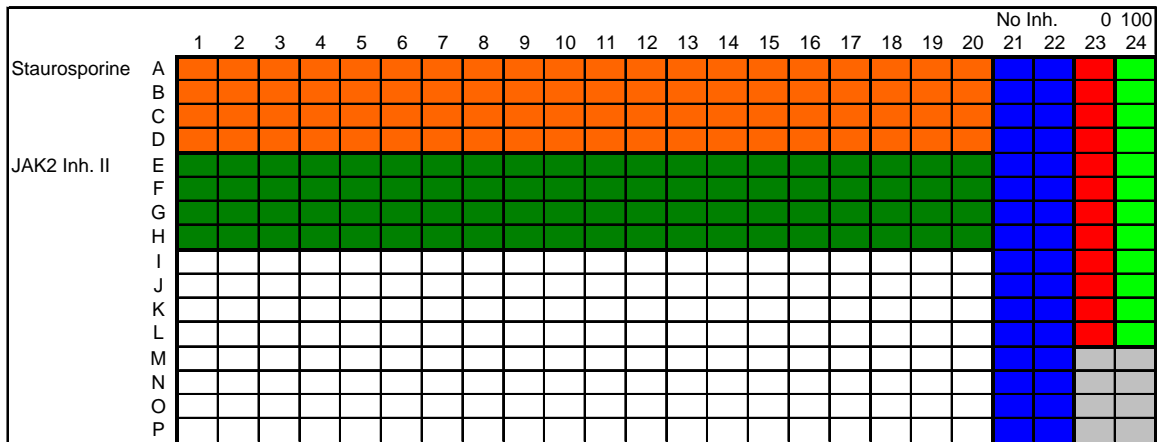
- from the next well to Column 2 Rows A through D and so on to set up an actual titration from 1  $\mu$ M final starting concentration of staurosporine).
8. Add 4  $\mu$ l Assay Buffer to Columns 21 and 22, as well as Columns 23 and 24, Rows A-L to bring all wells to a final volume of 40  $\mu$ l.
  9. Place TF-1 plate in incubator for 5 hours.

### Substrate Loading

10. Load cells as follows (note in both cases it is critical that cells are allowed to load at room temperature.):
  - For *Irf1-bla* TF-1 cells, 2.5 ml of 6X loading solution prepared by mixing 15  $\mu$ l CCF4-AM with 150  $\mu$ l Solution B, vortexing, then adding 2.335 ml of Solution C. Loading Solution mixed, and 8  $\mu$ l per well added to plate. Plate incubated at ROOM TEMP. for 2.5 hours.
  - For *Irf1-bla* HEL cells, 2.5 ml of 6X loading solution prepared by mixing 15  $\mu$ l CCF4-AM with 150  $\mu$ l Solution B, vortexing, and then adding 150  $\mu$ l Solution D and 2.185 ml Solution C. Loading Solution mixed, and 8  $\mu$ l per well added to plate. Plate incubated at ROOM TEMP. for 4 hours.

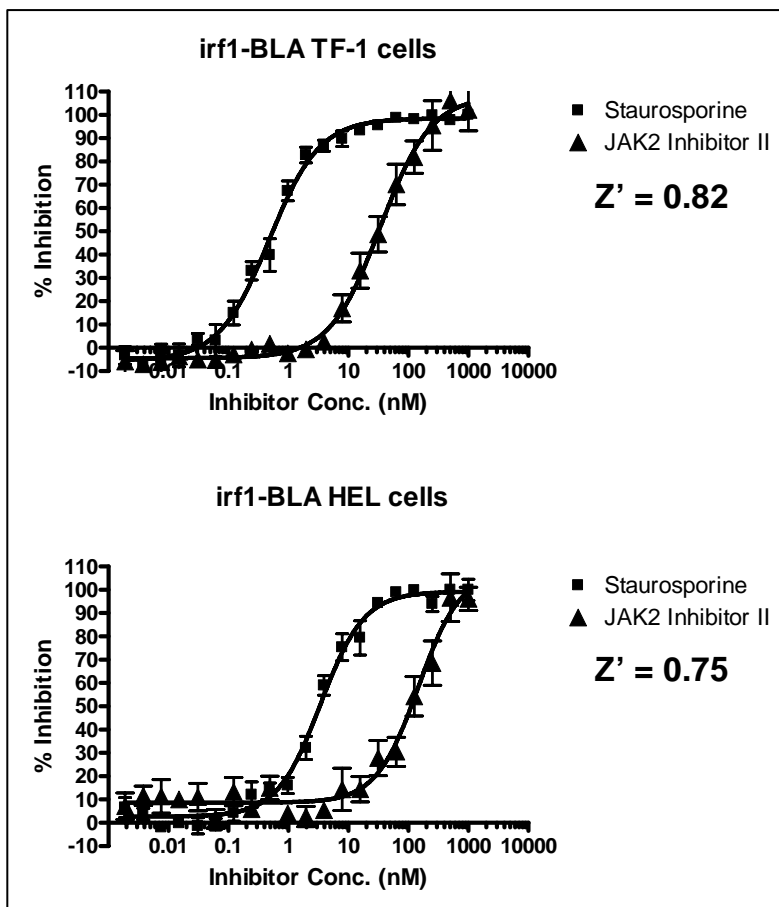
Setup Guide on the BMGLABTECH OPTIMA Microplate Reader

11. Read and analyze as directed in protocol.



**Figure 2: Assay Plate Schematic.** Compound titrations shown in Columns 1-20, Columns 21 and 22 prepared without any inhibitor as 0% inhibition controls, Column 23 prepared with no CM-CSF (unstimulated 0% activity control) for the TF-1 line, but the HEL line is constitutively active so Column 23 is untreated for HEL cells, and Column 24 is also untreated in both cases as a 100% activity control. Note 8 wells in gray in bottom right, which were prepared with out any inhibitor or cells, as controls for background subtraction.

D. Results



**Figure 1: GeneBLAzer® Assay.** GeneBLAzer® assay performed using the BMG LABTECH OPTIMA.