

**Modulus™ Microplate and Modulus™ II Microplate 96-well Data Conversion for Curve Fitting Data Analysis Software from Turner BioSystems**

Data generated using the Modulus™ Microplate and Modulus™ II Microplate instruments is in a format that is currently not compatible with using the **Load Raw Data** button that is one of the options for analyzing data using the Curve Fitting Data Analysis software provided by Turner Biosystems. If the **Load Raw Data** button is used a Null Reference Exception error message will appear on the screen as a result as can be seen in Figure #1.

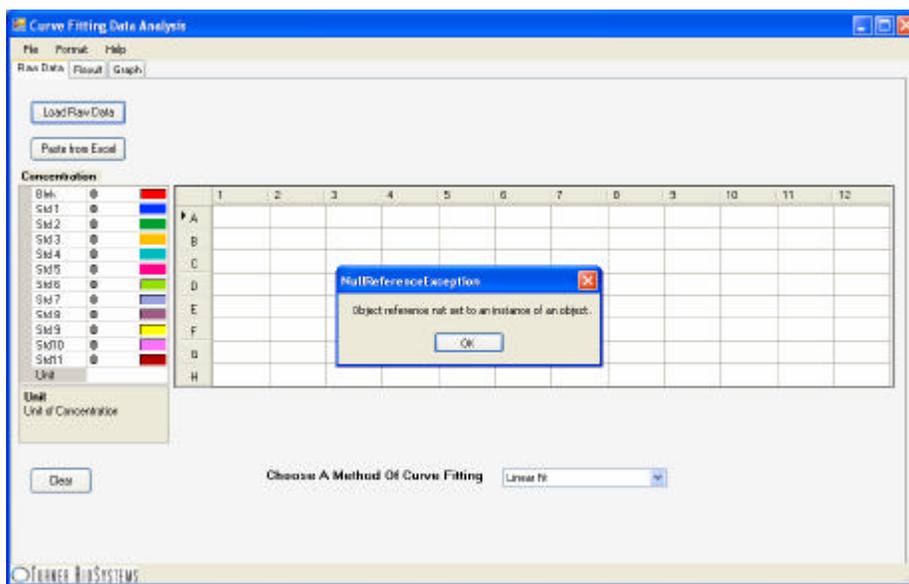


Figure #1: Null Reference Exception Error Message

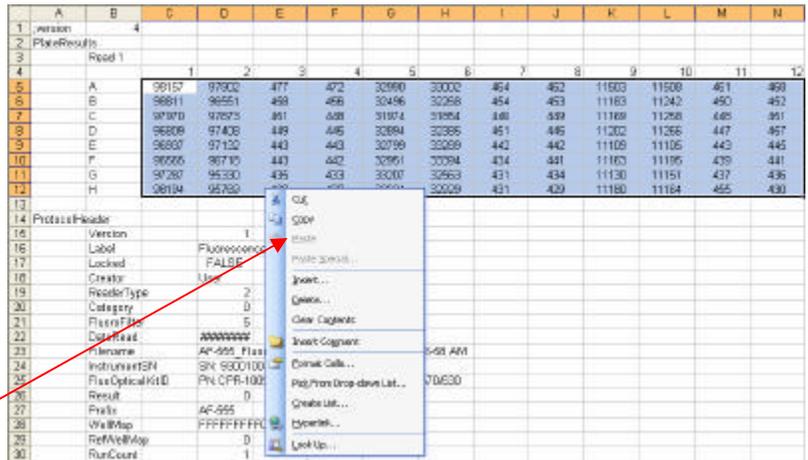
The following instructions are a simple exercise showing the method to convert 96-well data generated on the Modulus™ Microplate and Modulus™ II Microplate instruments into a format that is ideally suited for data plotting using the Curve fit data analysis software package from Turner BioSystems.

WRITTEN: J. Quast	DATE: April 15, 2009	Page 1 of 4		

Instructions for Pasting Data from Excel into Curve Fitting Data Analysis program

Step 1. Open .csv data file and highlight wells to be exported to Curve fitting software.

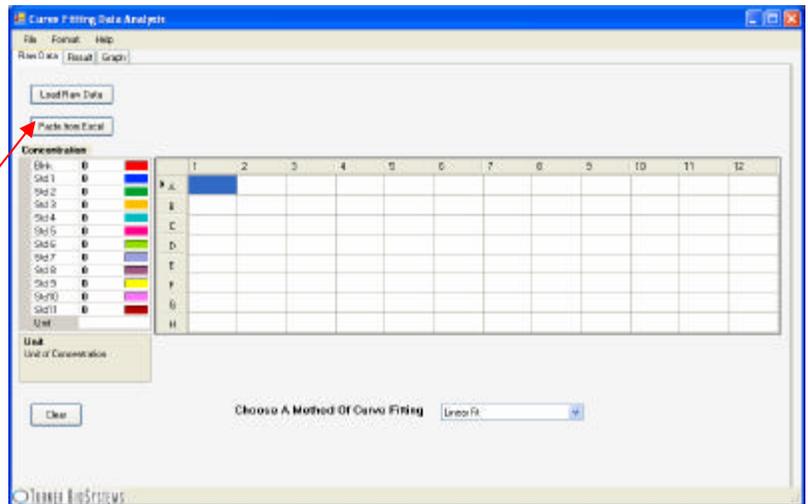
• **Note:** If wells were deselected during protocol set-up make certain that all X designations (deselected wells) and R designations (reference wells) are deleted before selecting data to be copied.



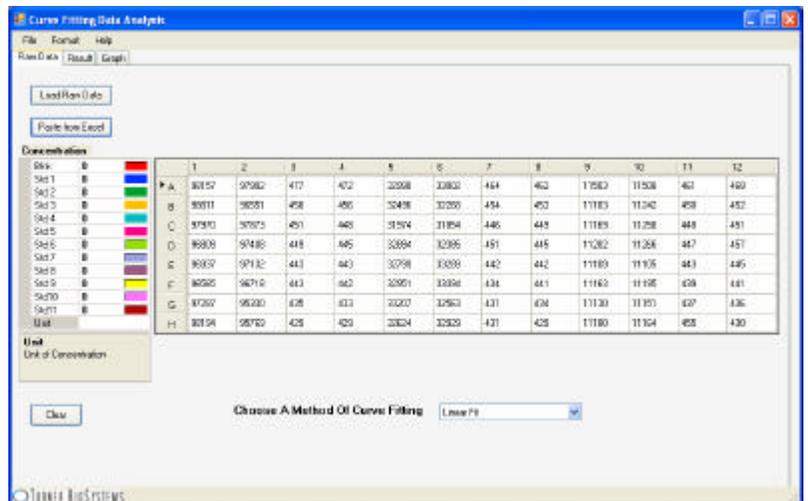
Step 2. Right-click on selected wells and select the Copy option from the pull down menu.

Step 3. Open the Curve Fitting Data Analysis program.

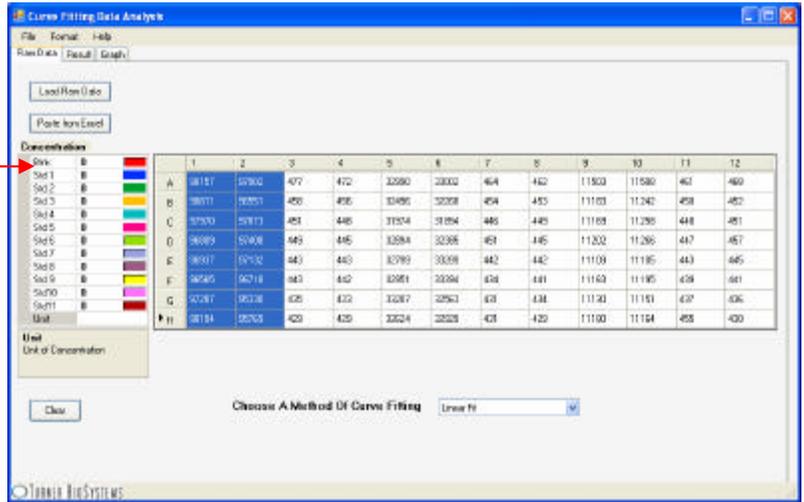
Step 4. Left click on well A1 and click on the **Paste from Excel** button to enter the raw data into the spreadsheet.



Step 5. With data Pasted into Curve Fitting Data Analysis spreadsheet, proceed with data reduction.



Step 6. Highlight the wells for Std 1 and click on the Std 1 cell to select. Repeat this for all other standards as well as the Blank (Blnk).



Step 7. Enter the values for the Standards in the cells to the right of the Std designations.

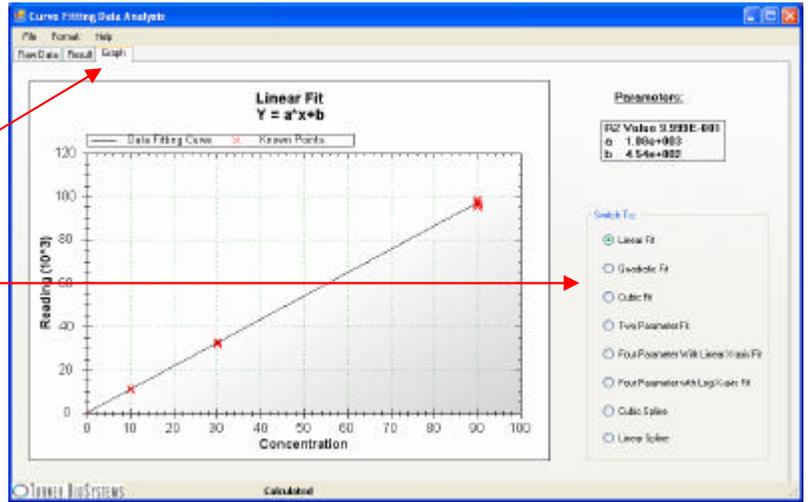


Step 8. To view the assay results, click on the **RESULT** tab.



Step 9. To view the assay results graph, click on the **GRAPH** tab.

Step 10. To view the assay results graph in other Curve fit parameters simply select from the box to the right of the graph.



**ABOUT TURNER BIOSYSTEMS, INC.**

Modulus is a trademark of Turner BioSystems. Orders for Turner BioSystems' products may be placed by:

Phone: (408) 636-2400  
 Toll Free: (888) 636-2401 or  
 Fax: (408) 737-7919

Contact us via our [contact form](#)  
 Internet: [www.turnerbiosystems.com](http://www.turnerbiosystems.com)

Mailing Address:  
 Turner BioSystems, Inc.  
 645 N. Mary Avenue  
 Sunnyvale, CA 94085

WRITTEN: J. Quast	DATE: April 15, 2009	Page 4 of 4		