

MetaXpress® 6.5 Software Guide

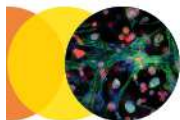
Plate Annotation

Date Revised 7/17/2019 Version B



What is plate annotation?

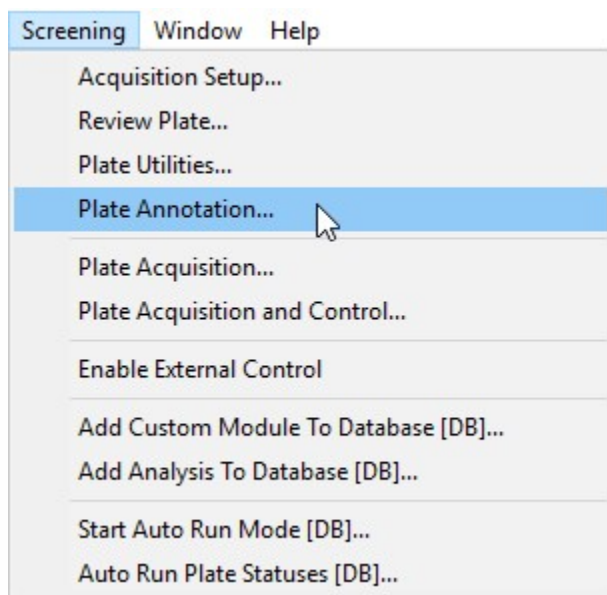
- Assigning sample information to specific wells of your plate:
 - Compound
 - Concentration
 - Category / group
- Annotation is specific for a plate
 - Applied across all measurement sets (analyses)
- Create annotation in MetaXpress/AcuityXpress, or import text file
- Batch annotation option for annotating multiple plates (for a screen)



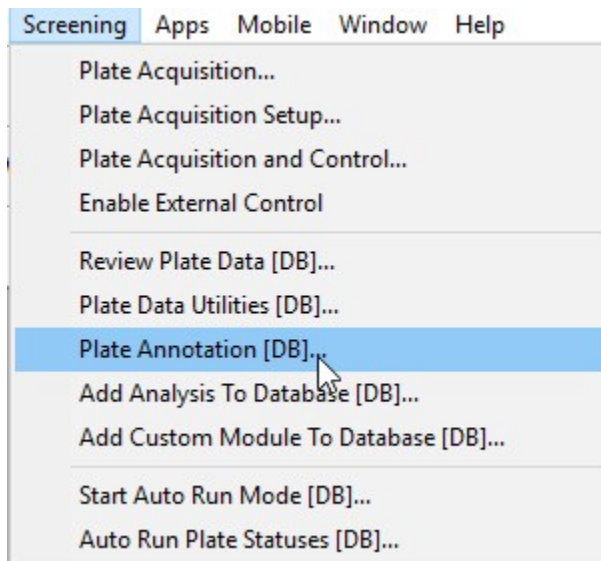
Editing Plate Annotation

- In MetaXpress version 6.5: **Screening > Plate Annotation**
- In the optional AcuityXpress software: **Manage > Plate Layouts**

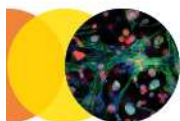
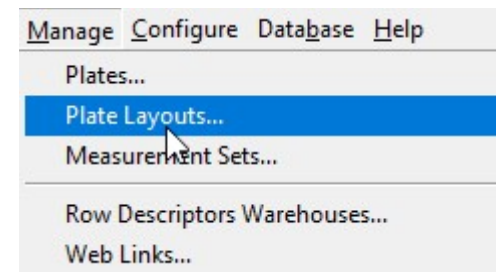
MX 6.5 Simplified Menu



MX 6.5 Standard Menu



AcuityXpress 2.1



Updating MX Simplified Menu

- If you do not see the **Plate Annotation** option in the simplified menu after updating to MetaXpress version 6.5, update the menu:

- Go to https://mdc.custhelp.com/app/answers/detail/a_id/20100 and download the latest version of the taskbar and documentation
- Go to **Control > Journal > Import Journal Suite** to import the taskbar, typically into C:\MX6\Taskbars
- Load the Main Taskbar
- Run the Taskbar Installer and select the option to **Configure simplified menu**
- Choose to **Install** the simplified menu.
- Select current group or all groups.
- Restart the MetaXpress program.

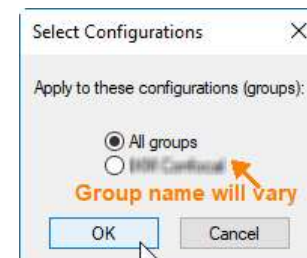
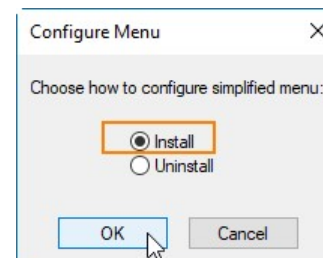
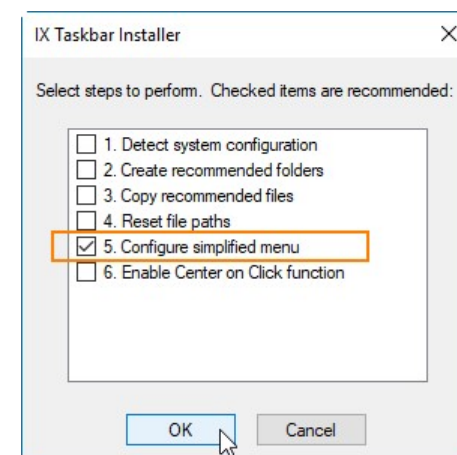
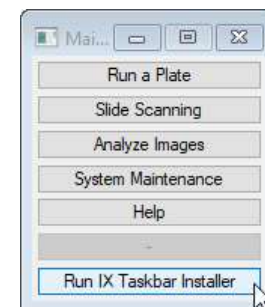


Plate Annotation interface

96 well (8 X 12), 1 compound, Untitled

Template: Untitled

Current Layout: Untitled

	01	02	03	04	05	06	07	08	09	10	11	12
A0	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M
B0	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M
C0	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M
D0	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M
E0	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M
F0	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M
G0	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M
H0	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M	Empty 0 μ M

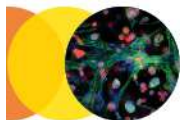
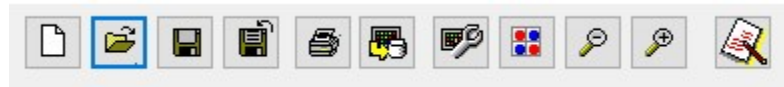
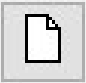









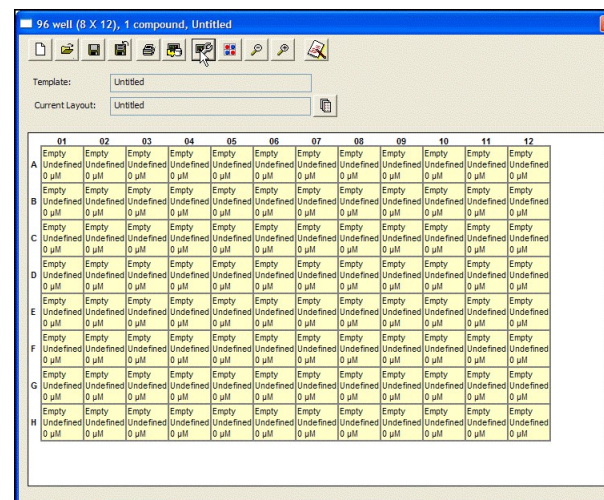


Plate Annotation options



-  • New (Ctrl + N)
-  • Open (Ctrl + O)
-  • Save (Ctrl + S)
-  • Save As (Ctrl + Shift + A)
-  • Print (Ctrl + P)
-  • Annotate Plate (Ctrl + D)
-  • Configure (Ctrl + Shift + C)
-  • Annotate Wells (Ctrl + Shift + S)
-  • Compact View
-  • Detailed View
-  • Batch Annotations (Ctrl + D)

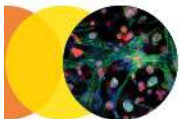


Typical workflows

Manually create plate layout

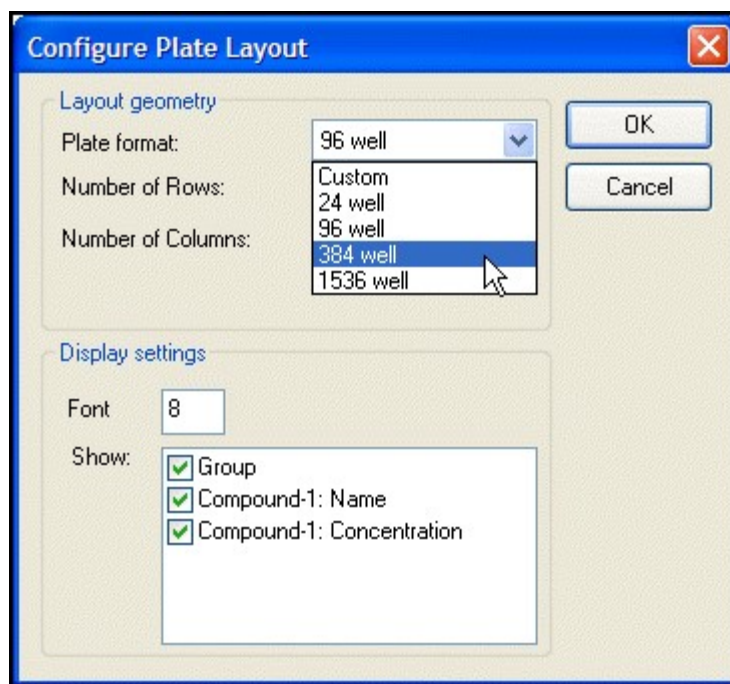


Import or open existing plate layout



Configure - Select plate format

- Select a pre-defined plate format or define a custom plate type.



Configure Plate Layout

Layout geometry

Plate format: 96 well

Number of Rows: Custom

Number of Columns: 384 well

OK

Cancel

Display settings

Font: 8

Show:

- Group
- Compound-1: Name
- Compound-1: Concentration



Annotate wells

- Highlight block of wells
- Click Annotate Wells icon
- Enter group, compound, and concentration information
- Repeat for other blocks of wells as necessary

	01	02	03	04	05	06	07	08	09	10	11	12
A	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM
B	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM
C	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM
D	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM
E	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM
F	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM
G	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM
H	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM



Example serial dilution

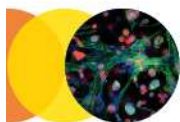
- Set up serial dilution from top to bottom
- All concentrations are in μM
- Top concentration is $50 \mu\text{M}$
- 5-fold dilutions down the rows
- Replicates in the columns

384 well (16 X 24), 1 compound, Untitled

Template: Untitled

Current Layout: Untitled

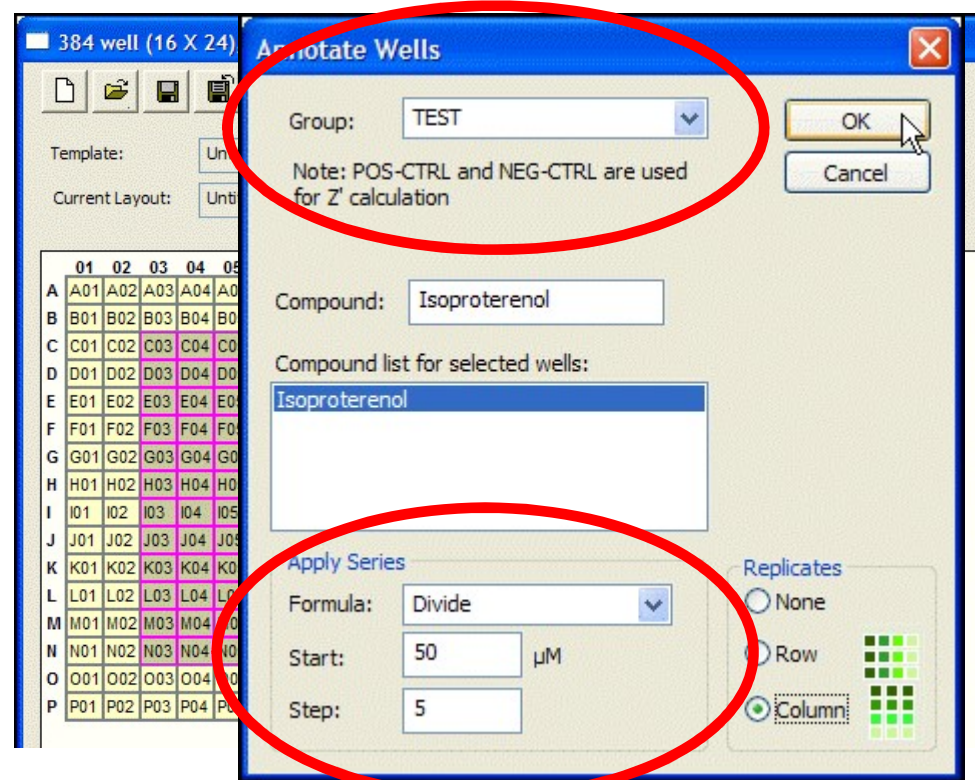
	01	02	03	04	05	06	07	08	09	10	11	12	13	14	15	16	17	18	19	20	21	22	23	24
A	A01	A02	A03	A04	A05	A06	A07	A08	A09	A10	A11	A12	A13	A14	A15	A16	A17	A18	A19	A20	A21	A22	A23	A24
B	B01	B02	B03	B04	B05	B06	B07	B08	B09	B10	B11	B12	B13	B14	B15	B16	B17	B18	B19	B20	B21	B22	B23	B24
C	C01	C02	C03	C04	C05	C06	C07	C08	C09	C10	C11	C12	C13	C14	C15	C16	C17	C18	C19	C20	C21	C22	C23	C24
D	D01	D02	D03	D04	D05	D06	D07	D08	D09	D10	D11	D12	D13	D14	D15	D16	D17	D18	D19	D20	D21	D22	D23	D24
E	E01	E02	E03	E04	E05	E06	E07	E08	E09	E10	E11	E12	E13	E14	E15	E16	E17	E18	E19	E20	E21	E22	E23	E24
F	F01	F02	F03	F04	F05	F06	F07	F08	F09	F10	F11	F12	F13	F14	F15	F16	F17	F18	F19	F20	F21	F22	F23	F24
G	G01	G02	G03	G04	G05	G06	G07	G08	G09	G10	G11	G12	G13	G14	G15	G16	G17	G18	G19	G20	G21	G22	G23	G24
H	H01	H02	H03	H04	H05	H06	H07	H08	H09	H10	H11	H12	H13	H14	H15	H16	H17	H18	H19	H20	H21	H22	H23	H24
I	I01	I02	I03	I04	I05	I06	I07	I08	I09	I10	I11	I12	I13	I14	I15	I16	I17	I18	I19	I20	I21	I22	I23	I24
J	J01	J02	J03	J04	J05	J06	J07	J08	J09	J10	J11	J12	J13	J14	J15	J16	J17	J18	J19	J20	J21	J22	J23	J24
K	K01	K02	K03	K04	K05	K06	K07	K08	K09	K10	K11	K12	K13	K14	K15	K16	K17	K18	K19	K20	K21	K22	K23	K24
L	L01	L02	L03	L04	L05	L06	L07	L08	L09	L10	L11	L12	L13	L14	L15	L16	L17	L18	L19	L20	L21	L22	L23	L24
M	M01	M02	M03	M04	M05	M06	M07	M08	M09	M10	M11	M12	M13	M14	M15	M16	M17	M18	M19	M20	M21	M22	M23	M24
N	N01	N02	N03	N04	N05	N06	N07	N08	N09	N10	N11	N12	N13	N14	N15	N16	N17	N18	N19	N20	N21	N22	N23	N24
O	O01	O02	O03	O04	O05	O06	O07	O08	O09	O10	O11	O12	O13	O14	O15	O16	O17	O18	O19	O20	O21	O22	O23	O24
P	P01	P02	P03	P04	P05	P06	P07	P08	P09	P10	P11	P12	P13	P14	P15	P16	P17	P18	P19	P20	P21	P22	P23	P24



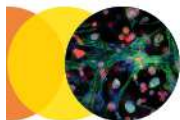
Example serial dilution

- Set up serial dilution from top to bottom
- All concentrations are in μM
- Top concentration is 50 μM
- 5-fold dilutions down the rows
- Replicates in the columns

Identify Controls or other Groups



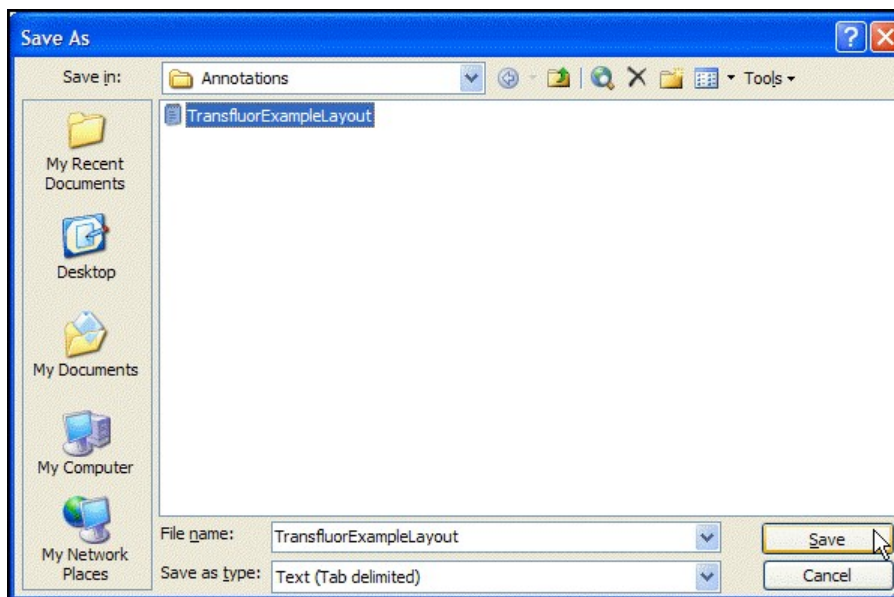
Serial Dilution



Or edit plate layout in Excel

- From Plate Annotation, select **Save As > File** to save an example plate layout file
- Edit the file in Excel
- Save the edited file as .txt (tab-delimited text)
- From Plate Annotation, select **Open > File** to import the annotations

	A	B	C	D	E
55	B14	Empty	Undefined	0	µM
56	B15	Empty	Undefined	0	µM
57	B16	Empty	Undefined	0	µM
58	B17	Empty	Undefined	0	µM
59	B18	Empty	Undefined	0	µM
60	B19	Empty	Undefined	0	µM
61	B20	Empty	Undefined	0	µM
62	B21	Empty	Undefined	0	µM
63	B22	Empty	Undefined	0	µM
64	B23	Empty	Undefined	0	µM
65	B24	Empty	Undefined	0	µM
66	C01	Empty	Undefined	0	µM
67	C02	Empty	Undefined	0	µM
68	C03	POS-CTRL	Isoproterer	50	µM
69	C04	POS-CTRL	Isoproterer	50	µM
70	C05	POS-CTRL	Isoproterer	50	µM
71	C06	TEST	TestCompi	50	µM
72	C07	TEST	TestCompi	50	µM
73	C08	TEST	TestCompi	50	µM
74	C09	Empty	Undefined	0	µM
75	C10	Empty	Undefined	0	µM



Select plate(s) to annotate

384 well (16 X 24), 1 compound, C:\Documents and Settings\paula\Desktop\2008-Imaging-CD\training\demodata\Annotations

Template: Untitled
Current Layout: Untitled

	01	02	03	04	05	06	07	08	09	10	11	12	
A	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	
B	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	
C	Empty Undefined 0 µM	Empty Undefined 0 µM	POS-CTRL Isoproterenol 50 µM	POS-CTRL Isoproterenol 50 µM	POS-CTRL Isoproterenol 50 µM	TEST Isoproterenol 10 µM	TEST Isoproterenol 10 µM	TEST Isoproterenol 10 µM	TEST Isoproterenol 10 µM	TEST Isoproterenol 10 µM	TEST Isoproterenol 2 µM	TEST Isoproterenol 2 µM	TEST Isoproterenol 2 µM
D	Empty Undefined 0 µM	Empty Undefined 0 µM	TEST Isoproterenol 0.4 µM	TEST Isoproterenol 0.4 µM	TEST Isoproterenol 0.4 µM	TEST Isoproterenol 0.08 µM	TEST Isoproterenol 0.08 µM	TEST Isoproterenol 0.08 µM	TEST Isoproterenol 0.08 µM	TEST Isoproterenol 0.016 µM	TEST Isoproterenol 0.016 µM	TEST Isoproterenol 0.016 µM	
E	Empty Undefined 0 µM	Empty Undefined 0 µM	TEST Isoproterenol 0.0032 µM	TEST Isoproterenol 0.0032 µM	TEST Isoproterenol 0.0032 µM	TEST TestCompound 0.0032 µM	TEST TestCompound 0.0032 µM	TEST TestCompound 0.0032 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	Empty Undefined 0 µM	
F	Empty Undefined 0 µM	Empty Undefined 0 µM	TEST Isoproterenol	TEST Isoproterenol	TEST Isoproterenol	TEST TestCompound	TEST TestCompound	TEST TestCompound	Empty Undefined	Empty Undefined	Empty Undefined	Empty Undefined	
G	Empty Undefined 0 µM	Empty Undefined 0 µM	TEST Isoproterenol	TEST Isoproterenol	TEST Isoproterenol	TEST TestCompound	TEST TestCompound	TEST TestCompound	Empty Undefined	Empty Undefined	Empty Undefined	Empty Undefined	
H	Empty Undefined 0 µM	Empty Undefined 0 µM	TEST Isoproterenol	TEST Isoproterenol	TEST Isoproterenol	TEST TestCompound	TEST TestCompound	TEST TestCompound	Empty Undefined	Empty Undefined	Empty Undefined	Empty Undefined	
I	Empty Undefined 0 µM	Empty Undefined 0 µM	TEST Isoproterenol	TEST Isoproterenol	TEST Isoproterenol	TEST TestCompound	TEST TestCompound	TEST TestCompound	Empty Undefined	Empty Undefined	Empty Undefined	Empty Undefined	
J	Empty Undefined 0 µM	Empty Undefined 0 µM	TEST Isoproterenol	TEST Isoproterenol	TEST Isoproterenol	TEST TestCompound	TEST TestCompound	TEST TestCompound	Empty Undefined	Empty Undefined	Empty Undefined	Empty Undefined	

Plate Dialog

- Plates
 - System Administrator [Creator Name - Plate Info]
 - 09/12/07 [Date Created - Plate Info]
 - 05/16/08 [Date Created - Plate Info]

Name [Plate Info]	Acquisiti...	Barcode ...	Creator ...	Date\Tim...
Demo dataset 5 Angiogenesis_PRIC...	Demo dat...	<NULL>	System A...	09/12/07...
dataset 1 adipogenesis_PRICKERT-U...	dataset 1...	<NULL>	System A...	09/12/07...
dataset 3 Neurite rinat_B_PRICKERT...	dataset 3...	18734J	System A...	09/12/07...
dataset 6 Translocation PKC_PRICK...	dataset 6...	<NULL>	System A...	09/12/07...
dataset 7 Apoptosis_PRICKERT-UCL...	dataset 7...	<NULL>	System A...	09/12/07...
TransfluorPitsBinned_PRICKERT-UCL...	Transflu...	<NULL>	System A...	09/12/07...
TransfluorVesidesBinned_PRICKERT...	Transflu...	<NULL>	System A...	09/12/07...

Plate Statistics

Plate Name: TransfluorVesidesBinned_PRICKERT-UCLT_29

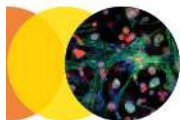
S: 72 W: 72 S: 72 C: 0 Cont...: 0 T: 4 M: 12

Select Cancel

Batch annotation

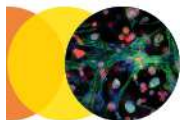
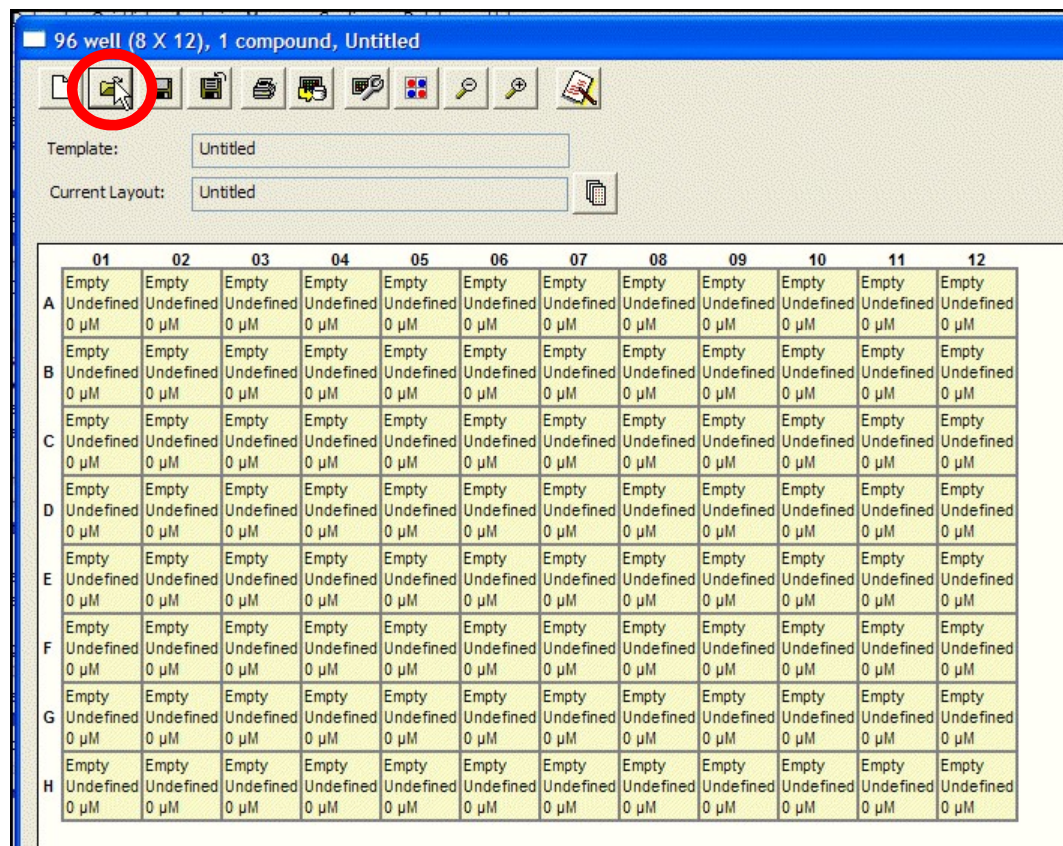
- Put multiple plates layouts in one text file
- Identify individual plates through one of the following:
 - Barcode
 - Plate Name
 - Plate ID
- Set “ID_Name” and “ID_Value” accordingly

\$HeaderStart\$			
TemplateName=Untitled			
TemplateDesc=			
\$HeaderEnd\$			
\$LayoutHeaderStart\$			
Rows=16			
Columns=24			
Format=0			
Compounds=1			
Description=ND0011637			
Name=Transfluor LOPAC Agonist ND0011637			
ID_Name=Barcode			
ID_DisplayName=			
ID_Value=ND0011637			
\$LayoutHeaderEnd\$			
\$LayoutDataStart\$			
WELL	GROUP	COMPOUND	CONCENT UNIT1
A01	Empty	Undefined	0 µM
A02	Empty	Undefined	0 µM
A03	TEST	DL-alpha-N	10 µM
A04	TEST	N-Acetyl-L	10 µM
A05	TEST	6-Methoxy	10 µM
A06	TEST	6-Aminohe	10 µM
A07	TEST	Acetamide	10 µM
A08	TEST	Altretamin	10 µM
A09	TEST	Amantadin	10 µM



Batch Annotation

- Open text file in the Plate Layout Manager

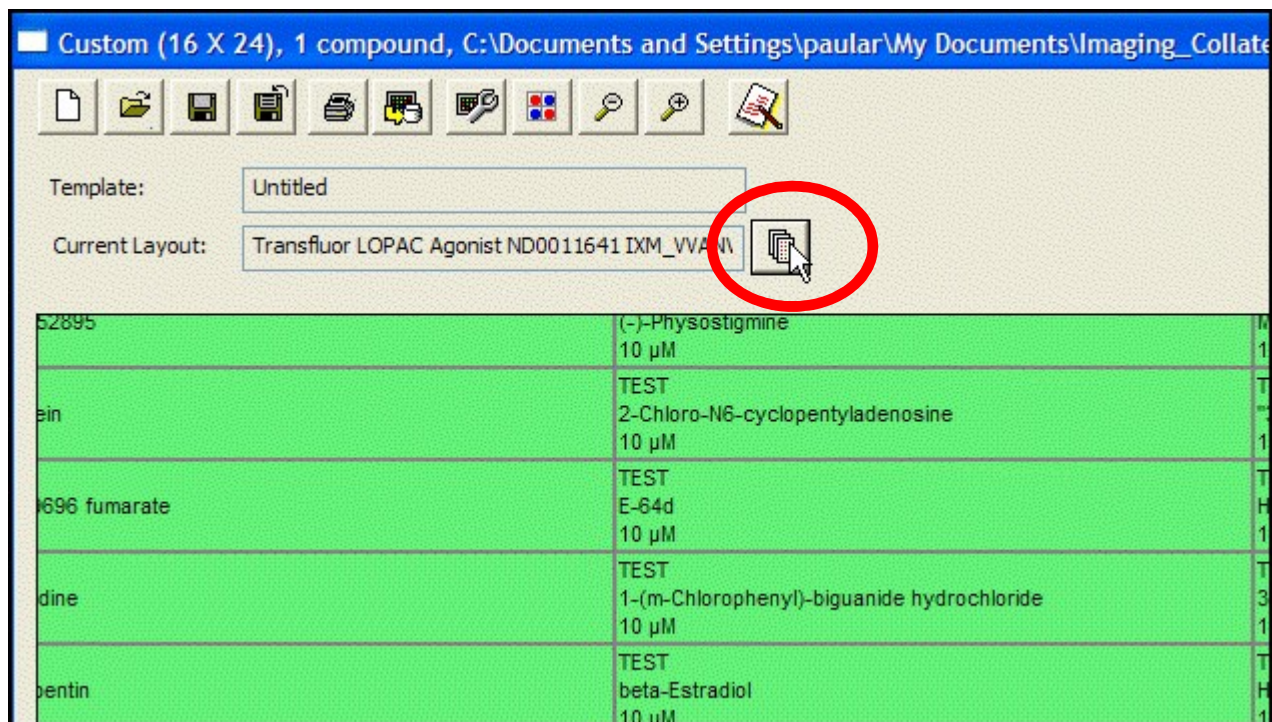


Batch Annotation – opening layout

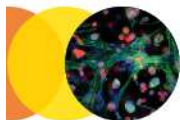
The image shows a software interface with three overlapping windows:

- Warning Dialog:** A dialog box with a red 'X' icon and the text "Do you want to save changes made to the plate layout?". It has three buttons: "Yes", "No", and "Cancel".
- Open Layout Dialog:** A dialog box titled "Open Layout" with the text "Open layout from". It has four radio button options: "File" (selected), "Template in the database", "Plate in the database", and "Measurement Sets associated with a plate (Use this option if plate annotations are stored in the measurement sets that were imported from file or annotated with previous version of the software)". It has "OK" and "Cancel" buttons.
- File Explorer:** A window showing a folder named "annotations". The file list includes:
 - Transflur_Agonist_A_Layout.txt
 - Transflur_Agonist_B_Layout.txt
 - Transflur_Agonist_C_Layout.txt
 - Transflur_Agonist_D_Layout.txt
 - Transflur_Agonist_Screen_b.txt
 - Transflur_Agonist_Screen_Batch.txt (highlighted)
 - Transflur_Agonist_Screen_ND0011637.txt
 - Transflur_Agonist_Screen_ND0011641.txt
 - Transflur_Agonist_Screen_ND0011645.txt
 - Transflur_Agonist_Screen_ND0011649.txtThe "File name" field contains "Transflur_Agonist_Screen_Batch.txt" and the "Files of type" field is set to "Plate Layout files (*.txt)". "Open" and "Cancel" buttons are visible.

Batch annotation



- Click Select icon to view different layouts in batch template



Batch annotation

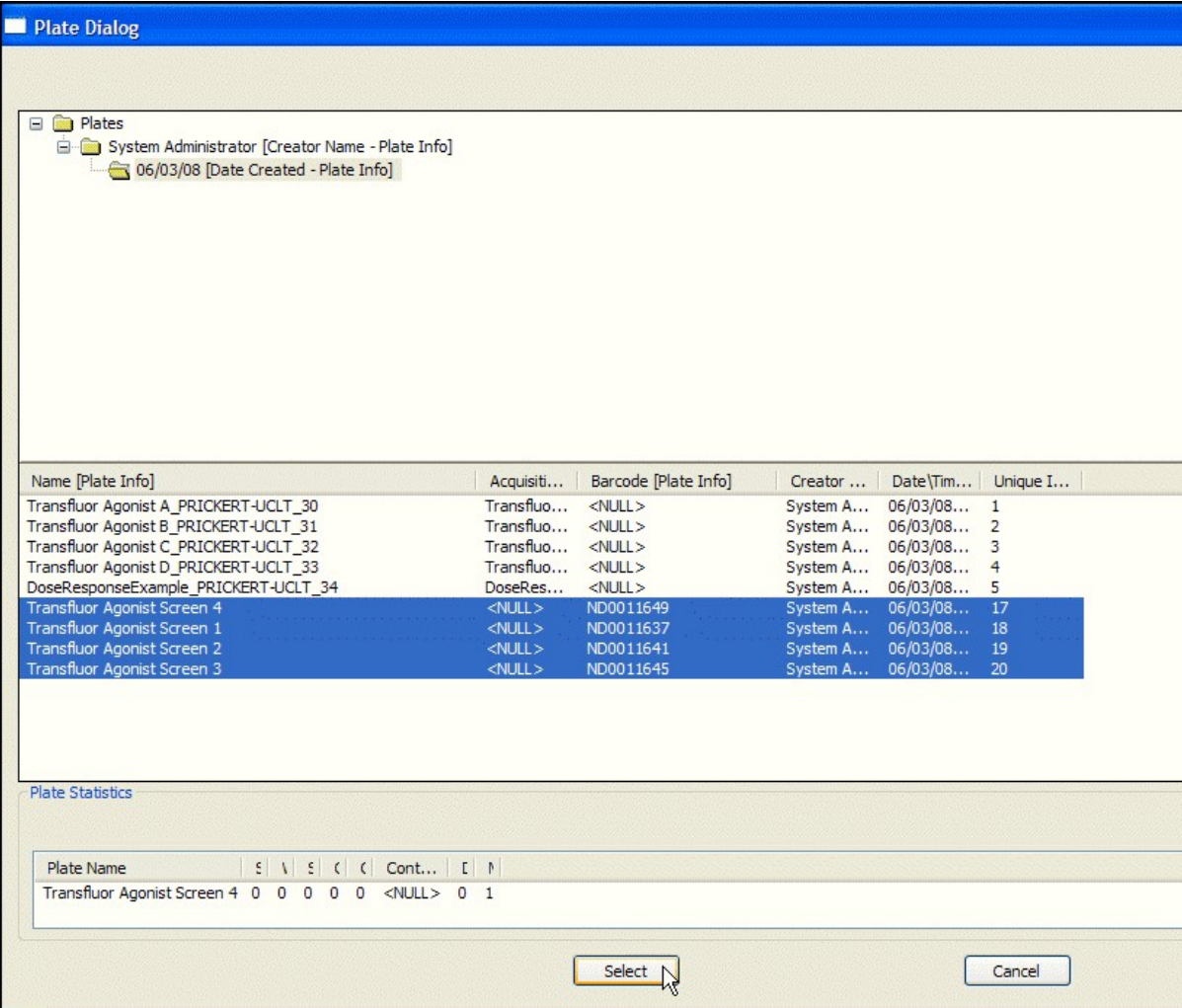
- Use the **Batch Annotations** button

Compound	Test	Control/Compound
Amiloride dicyanide	Iodoacetamide 10 µM	Isoproterenol 1 µM
	TEST	POS-CTRL
	Etoposide 10 µM	Isoproterenol 1 µM
	TEST	POS-CTRL
	HA-100 10 µM	Isoproterenol 1 µM
	TEST	POS-CTRL
Terbutaline-N-oxide*	ET-18-OCH3 10 µM	Isoproterenol 1 µM
	TEST	POS-CTRL
	Ipratropium bromide 10 µM	Isoproterenol 1 µM
	TEST	POS-CTRL
Terbutaline*	Etazolate hydrochloride 10 µM	Isoproterenol 1 µM
	TEST	POS-CTRL
	Idarubicin 10 µM	Isoproterenol 1 µM



Batch annotation

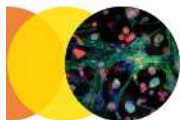
- Select all plates for batch annotation
- Use Ctrl or Shift to multi-select



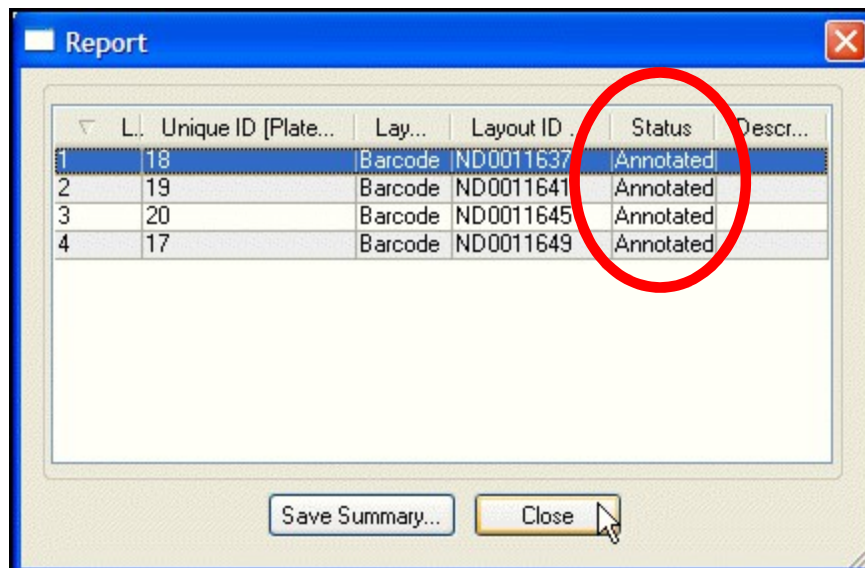
The screenshot shows the 'Plate Dialog' window. The top section is a tree view showing a folder structure: 'Plates' > 'System Administrator [Creator Name - Plate Info]' > '06/03/08 [Date Created - Plate Info]'. Below this is a table with the following columns: Name [Plate Info], Acquisiti..., Barcode [Plate Info], Creator ..., Date\Tim..., and Unique I... The table contains 10 rows of data, with the last five rows highlighted in blue. Below the table is a 'Plate Statistics' section with a table showing the statistics for the selected plate.

Name [Plate Info]	Acquisiti...	Barcode [Plate Info]	Creator ...	Date\Tim...	Unique I...
Transfluor Agonist A_PRICKERT-UCLT_30	Transflu...	<NULL>	System A...	06/03/08...	1
Transfluor Agonist B_PRICKERT-UCLT_31	Transflu...	<NULL>	System A...	06/03/08...	2
Transfluor Agonist C_PRICKERT-UCLT_32	Transflu...	<NULL>	System A...	06/03/08...	3
Transfluor Agonist D_PRICKERT-UCLT_33	Transflu...	<NULL>	System A...	06/03/08...	4
DoseResponseExample_PRICKERT-UCLT_34	DoseRes...	<NULL>	System A...	06/03/08...	5
Transfluor Agonist Screen 4	<NULL>	ND0011649	System A...	06/03/08...	17
Transfluor Agonist Screen 1	<NULL>	ND0011637	System A...	06/03/08...	18
Transfluor Agonist Screen 2	<NULL>	ND0011641	System A...	06/03/08...	19
Transfluor Agonist Screen 3	<NULL>	ND0011645	System A...	06/03/08...	20

Plate Name	ε	\	ε	<	<	Cont...	Γ	Π
Transfluor Agonist Screen 4	0	0	0	0	0	<NULL>	0	1



Batch annotation



L.	Unique ID [Plate...]	Lay...	Layout ID	Status	Descr...
1	18	Barcode	ND0011637	Annotated	
2	19	Barcode	ND0011641	Annotated	
3	20	Barcode	ND0011645	Annotated	
4	17	Barcode	ND0011649	Annotated	

- Check the resulting report to verify each plate is “Annotated”



Updating plate annotations

- You can update the plate annotation at any point, except:
 - If you have made a dataset from that plate in the optional AcuityXpress software
 - In this case, changes to the plate annotation would invalidate analysis results
- You can update plate annotation after viewing curve fit or peak analysis results in the MetaXpress software

