

MetaXpress® 6 Software Guide

Running an Analysis in Review Plate Data



Date Revised 07/28/15 Version B

2012-2015. Trademarks property of Molecular Devices, LLC or their respective owners. For research use only. Not for use in diagnostic procedures.

Chapter Purpose

The purpose of this chapter is to guide the user through running an analysis in the **Review Plata Data** dialog for single plane, Timelapse, or Z Series data sets.

The **Review Plate Data** dialog allows you to run an analysis on all wells, selected wells, selected site, and/or selected Time points or Z Steps on the plate displayed.

This chapter <u>will not</u> go over how to set up an analysis or describe batch analysis. Refer to corresponding chapters for details.





Running an Analysis in Review Plate Data

- 1. Open Review Plata Data
 - In the main toolbar click on



OR

- Under the Screening menu, select Review Plate Data
- 2. Click on the Select Plate button
- 3. Browse through the folders to open the plate of interest

Vavelengths:	Data v	iew:	: N	/ell a	arra	nge	mer	nt			•										F	nint	Tal	ble	
DAPI		01	02	03	04	05	06	07	08	09	10	11	12	13	14	15	16	17	18	19	20	21	22	23	24
FITC	A																								
	В																								
	С			•																					
	D			•																					
	E			•																					
	F			•														_							
	G			•																					
	н			•																					
				•																					
	J			•																					
Legend	K			•																					
Not acquired	L			•				_																	
Acquired, not measured	м			•																					
Displayed well	N			·																					
Part of montage	0							_																	
Selected wells	Р																								

Select Plate for Review	Plate Info]			
			I	1
Plates System Administrator [Creator Name - O4/20/15 [Date Created - Plate Info] O5/21/15 [Date Created - Plate Info] O6/04/15 [Date Created - Plate Info]	Plate Info]			
Name [Plate Info]	Acquisiti	D 1		
	Acquisiti	Barcode	Creator	Date\Ti
EX1NuclearFoci_AMSNVL-C0H8KV1_1	EX1Nucl	<null></null>	Creator System	Date\Ti 04/20/15
EX1NuclearFoci_AMSNVL-C0H8KV1_1 EX2TransfluorVesicles_AMSNVL-C0H8KV1_2	EX1Nucl EX2Tran	<null> <null></null></null>	System System	Date\Ti 04/20/15 04/20/15
EX1NuclearFoci_AMSNVL-C0H8KV1_1 EX2TransfluorVesicles_AMSNVL-C0H8KV1_2 EX3ThreeCHCellScoring_AMSNVL-C0H8KV1_3	EX1Nucl EX2Tran EX3Thre	<pre>And And And And And And And And And And</pre>	System System System	Date\Ti 04/20/15 04/20/15 04/20/15





Run an Analysis

- 4. Select the Run Analysis tab in the Review Plate Data dialog
- 5. Select the appropriate **Analysis** and **Settings** from the drop-down menus
 - Application modules will appear in <> brackets
 - Journals and Custom Modules do not have <> brackets
 NOTE If the selected analysis is a custom module, the Configure Settings button will revert Configure
 Custom Module. If the select analysis is a journal, the button will revert to Run Setup for Analysis.
- 6. Make sure Log into the database is enabled

	Montage: 1 🚔 x 3 荣]	
Display R	un Analysis Measurements Graph			1	
Analysis:	<transfluor></transfluor>	•	Configure Settings		
Settings:	Transfluor Vesicles	Edit List	Create Custom Module		After analysis has been
Setting	Transfluor Vesciles 05062015		A		completed on the first well,
description:			Ŧ		and approximate time to
	Remaining Time:		Run on all wells		finish will appear.
	0.6 min		Run on selection		
	Cog int	to the database	Run on displayed site		





Run an Analysis on All Wells – Single Plane

To run the select analysis on the all the images in the plate:

- 7. Click on the **Run on all wells** button
 - If the plate contains sites, analysis will be ran on all sites in all wells

	Montage: 1 x 3					
Display R	un Analysis Measurements Graph					
Analysis:	<transfluor></transfluor>	•	Configure Settings			
Settings:	Transfluor Vesicles 🔹	Edit List	Create Custom Module			
Setting description:	Transfluor Vesciles 05062015		* *			
	Remaining Time:	Remaining Time:				
	0.6 min	0.6 min				
	Log into	the database	Run on displayed site			





Run an Analysis on Selected Wells – Single Plane

To run an analysis on selected wells

7. In the plate view, right-click on wells of interest to highlight them in green

NOTE If an analysis has previously been ran on this plate, deselect the Heat Map option on the Measurements tab to easily see selected wells

- 8. Click on the **Run on selection** button
 - If the plate contains multiple sites, analysis will be ran on all sites in selected wells







Run an Analysis on Displayed Site – Single Plane

To run an analysis on a single site (displayed image)

7. In the thumbnail montage, click on well of interest. A red border will appear around the selected well in the plate grid.

NOTE DO NOT click in the plate grid as this will only change the wells displayed in the thumbnail montage

- 8. Click on the **Run on displayed site** button
 - If the plate contains multiple sites, analysis will be ran on the site highlighted in gray in the site section







Run an Analysis on Timelapse Data

If the image data set contains Timelapse images:

- 7. Select the appropriate selection under the **Time points** section:
 - All time points: all time points in the data set will be analyzed
 - **Time point range**: only the specified contiguous time points in the spin boxes will be analyzed
 - Selected time point: only the time point selected in the Time points spin box below the plate grid will be analyzed
 - **Stack of all time points**: this option is only available when a journal is selected for analysis

Display F	Montage: 2 Run Analysis Measurements Gra	★x 1 ★ Time points: 2 ★ of 6	
Analysis:	<transfluor></transfluor>	•	Configure Settings
Settings:	Transfluor Vesicles	▼ Edit List	Create Custom Module
Setting description	Transfluor Vesciles 05062015		* *
Time poir All tim Time p Select Stack	nts: le points point range 1 1 6 0 ted time point c of all time points	☑ Log into the database	Run on all wells Run on selection Run on displayed site





Run an Analysis on Timelapse Data

- 8. As with single plane data sets, choose the appropriate selection for running the analysis:
 - Run on all wells: analysis will be ran on all sites/wells
 - **Run on selection**: analysis will be run on all selected sites/wells. Right click on wells of interest in the plate grid to highlight in green.
 - Run on displayed site: analysis will be run on the site highlighted in the plate grid with a red border

NOTE For more details, refer to sections on running an analysis on a single plane.

	Montage: 2 x 1	Time points: 2 🌩 of 6	
Display R	un Analysis Measurements Graph		
Analysis:	<transfluor></transfluor>	•	Configure Settings
Settings:	Transfluor Vesicles	▼ Edit List [Create Custom Module
Setting description:	Transfluor Vesciles 05062015		A
Time point All time Time p Selecter Stack	ts: points oint range 1 1 6 1		Run on all wells Run on selection
U Stack	or all time points	Log into the database	Run on displayed site





Run an Analysis on Z Series Data

If the image data set contains Z Series images:

- 7. Select the appropriate selection under the **Z steps** section:
 - All Z steps: all Z steps in the data set will be analyzed
 - **Z step range**: only the specified contiguous Z steps range in the spin boxes will be analyzed
 - Selected Z step: only the Z step selected in the Z steps spin box below the plate grid will be analyzed
 - Stack of all Z steps: this option is only available when a journal is selected for analysis
 - **2D Projection**: only the 2D projection image in the data set will be analyzed

	Montage: 2	⇒ x 1 🜩	Z steps:	2	🖨 of 5		2D projection		•
Display F	Run Analysis Measurements Gra	ph							
Analysis:	<transfluor></transfluor>					•	Configure Se	ttings	
Settings:	Transfluor Vesicles			•	Edit List	. [Create Custon	n Module	
Setting description:	Transfluor Vesciles 05062015							^ ~	
Z steps: All Z steps:	teps						Run on a	ll wells	וו
Select	ed Z step						Run on se	election]
© 2D pro	or all ∠ steps		🔽 L	og into	the databa	ase	Run on disp	layed site]





Run an Analysis on Z Series Data

- 8. As with single plane data sets, choose the appropriate selection for running the analysis:
 - Run on all wells: analysis will be ran on all sites/wells
 - **Run on selection**: analysis will be run on all selected sites/wells. Right click on wells of interest in the plate grid to highlight in green.
 - Run on displayed site: analysis will be run on the site highlighted in the plate grid with a red border

NOTE For more details, refer to sections on running an analysis on a single plane.

	Montage: 2 x 1	Z steps: 2 🔹 of 5	2D projection	
Display R	lun Analysis Measurements Graph			
Analysis:	<transfluor></transfluor>		Configure Settings	
Settings:	Transfluor Vesicles	▼ Edit L	ist Create Custom Module	
Setting description:	Transfluor Vesciles 05062015			^ -
Z steps: All Z st Z step Selecte	teps range 1 ∳ 5 ∳ ed Z step		Run on all wells Run on selection	
 Stack 2D proj 	of all Z steps jection	🗹 Log into the da	tabase Run on displayed sit	e





Support Resources

- F1 / HELP within MetaXpress® Software
- Support and Knowledge Base: <u>http://mdc.custhelp.com/</u>
- User Forum: http://metamorph.moleculardevices.com/forum/
- Request Support: <u>http://mdc.custhelp.com/app/ask</u>
- Technical Support can also be reached by telephone:
 - 1 (800) 635-5577
 - Select options for Tech Support → Cellular Imaging Products → ImageXpress Instruments





MOLECULAR DEVICES

ADVANCING PROTEIN AND CELL BIOLOGY

For research use only. Not for use in diagnostic procedures.