



MetaXpress[®] 6 Software Guide

Enabling Adaptive Acquisition[™]

Date Revised 06/08/15 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 enabling the **Adaptive Acquisition** feature in **Plate Acquisition Setup**.

When the **Adaptive Acquisition** feature is enabled, the system images the minimum number of sites to reach the user-defined target number of cells within each well. It is particularly useful for assays in which the cell seeding or distribution is uneven from well to well.



Adaptive Acquisition Overview

When to use

- You need a minimum number of cells for statistical significance, but there is uneven cell seeding from well to well
 - Translocation (% positive measurement)
 - Cell cycle (% G1, % M, etc.)
- To detect rare events
- Rare GFP-positive cells (any present)

How it works

- Specify a minimum number of cells to be counted per well
- System acquires sites and counts cells until minimum cell count is reached
- System moves stage to the next well

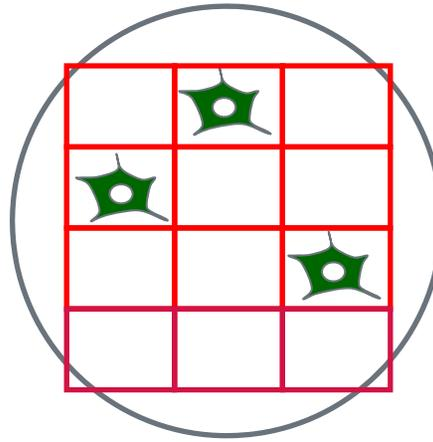


How Does Adaptive Acquisition Work?

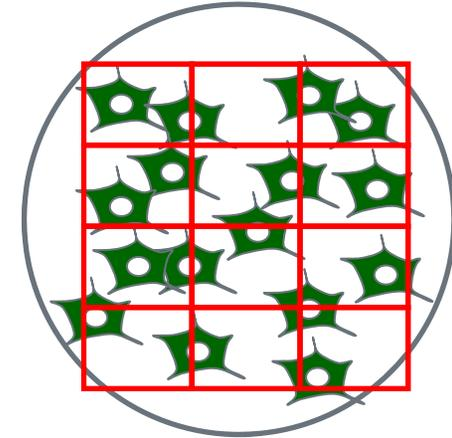
Without Adaptive Acquisition

- Target # of cells: 3
- Always acquire all 12 sites

12 Sites Acquired



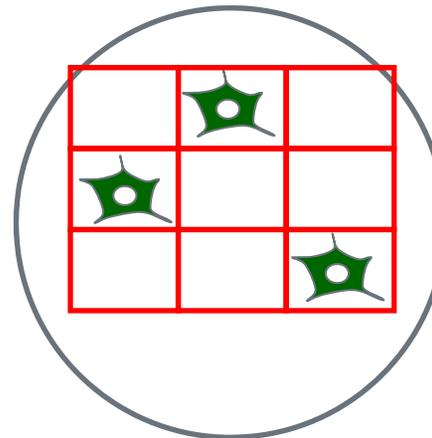
12 Sites Acquired



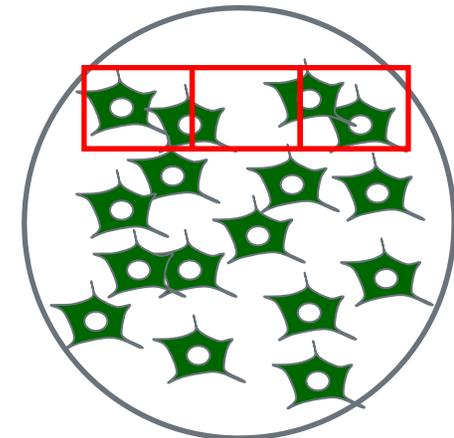
With Adaptive Acquisition

- Target # of cells: 3
- Acquire variable number of sites to reach target cell count

9 Sites Acquired



3 Sites Acquired



Enabling Adaptive Acquisition

1. Open Plate Acquisition Setup

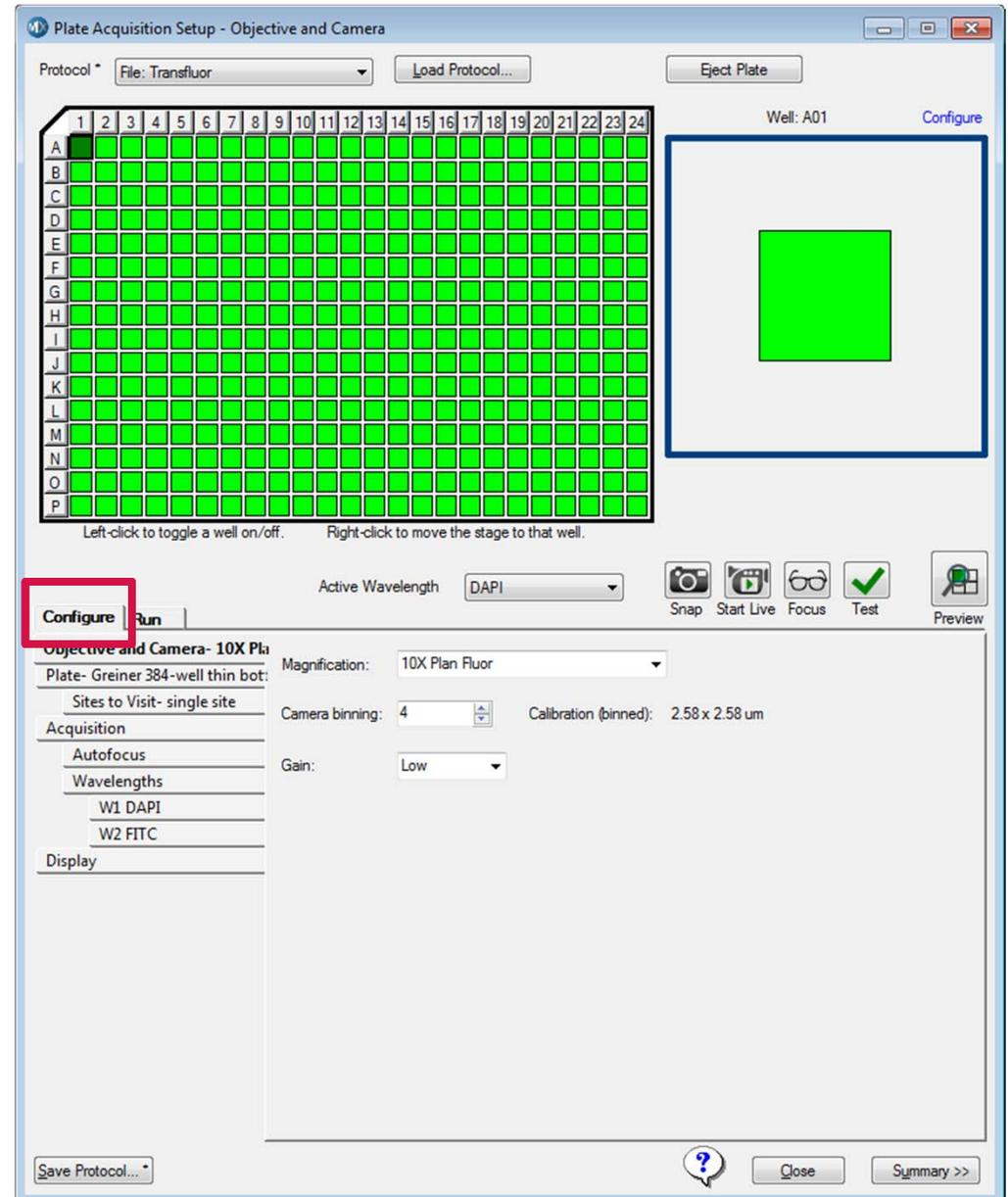
- In the main toolbar click on



OR

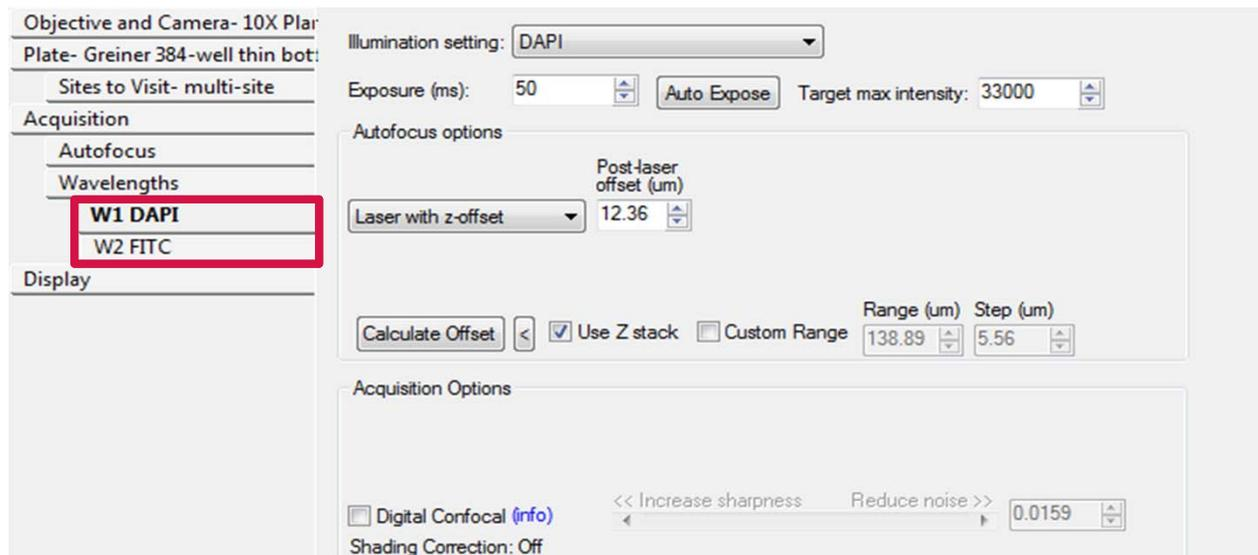
- Under the **Screening** menu, select **Plate Acquisition Setup**

2. Select the **Configure** tab



Enabling Adaptive Acquisition

3. Select the appropriate **W** tab (wavelength) for which you will use to count cells
4. Configure the wavelength settings as you would for acquisition
 - i. Select correct **Illumination Setting** from the drop-down menu
 - ii. Calculate **Focus offset**
 - iii. Determine **Exposure time**
5. After you have optimized settings, right click on a site/well with representative cells as this this position will be used to test settings



Enabling Adaptive Acquisition

5. Select the **Site to Visit** tab
6. Under the **Site Options** section, select **Adaptive Acquisition**
 - The Adaptive Acquisition section appears at the bottom of the dialog box
7. Select the appropriate (max) **number of sites** and **spacing**
**NOTE* The system will acquire at most the number of sites indicated in this section even if the cell count per well has not been reached*

Objective and Camera- 40X S PI
Plate- Greiner 384-well thin bot:

Sites to Visit- adaptive

Acquisition

Autofocus

Wavelengths

W1 DAPI

W2 FITC

Display

Site Options

Single site

Fixed number of sites

Adaptive acquisition

Multi-well

Custom field of view (%):
X: 50 Y: 50

Well size: 11 mm²
Number of sites: 81
90.18% Well Coverage

Site/image size: 348.19 x 348.19 μm

Acquires sites based on the number of cells per well

Columns: 9 Spacing (μm) 0

Rows: 9 Spacing (μm) 0

Tile sites

Fit sites to well

Overlap sites 10%

Adaptive Acquisition

Minimum sites to visit: 2

Test Segmentation

Wavelength: W1 - DAPI

Nuclei count: 0

Approximate width: 5 to 10 μm

Intensity above local background: 100 gray levels

Cell Count per well: 50



Configuring Adaptive Acquisition Settings

8. Enter the **Minimum number of sites to visit**: system will acquire at least this number of sites per well even if the target cell count has already been reached
9. Select the cell counting **Wavelength**: should be the first wavelength acquired and is often a nuclear marker
**NOTE*: This must be the same wavelength chosen on the W1 tab*
9. Enter the cell counting parameters (see next section for details):
 - **Approximate minimum and maximum width** of cells
 - **Intensity above local background** for the cell marker
 - Specify target **Cell Count per well**: the minimum number of cells desired by the user

Adaptive Acquisition

Minimum sites to visit: 2

Test Segmentation

Nuclei count: 0

Wavelength: W1 - DAPI

Approximate width: 5 to 10 μm

Intensity above local background: 100 gray levels

Cell Count per well: 50



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

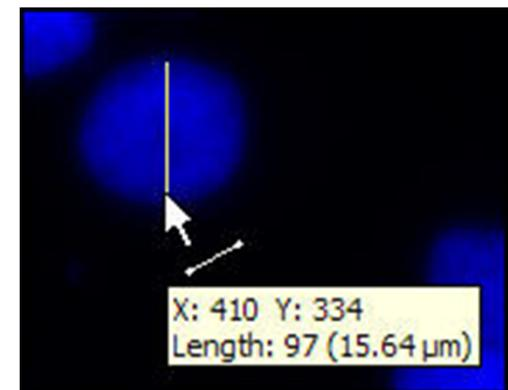
Determining Approximate Width Settings

Set the **Approximate min width** and **Approximate max width** for the range of cells (or nuclei) that you want to detect.

- Select the line region from the **Region Tools** toolbar
- Click and drag across a representative small and large cell; a tooltip will show the length of the line.

NOTE Do not click the image again as this will cause the tooltip to disappear. If the tooltip disappears, repeat the drawing procedure as the lines will not affect any image settings.

- The width is the short axis of a cell (in μm)
- Much smaller or much larger cells will be ignored



Effects of Varying Width Settings

Molecular Devices recommends only changing one parameter at a time in order to determine optimal settings.

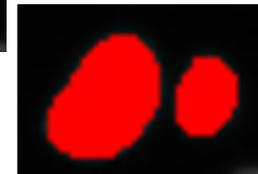
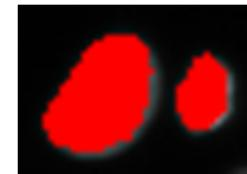
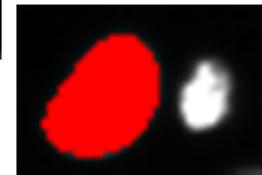
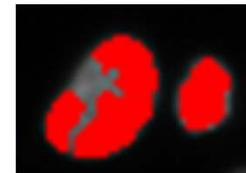
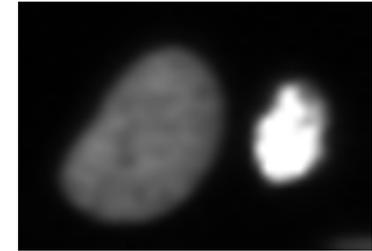
Min width too small: splits nuclei

Min width too large: omits smaller nuclei

Max width too small: may shrink nuclear boundaries

Max width too large: may slightly enlarge nuclear boundaries

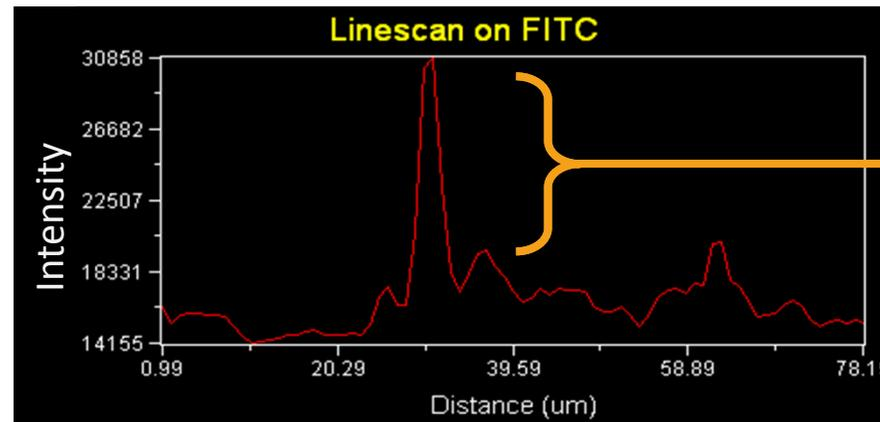
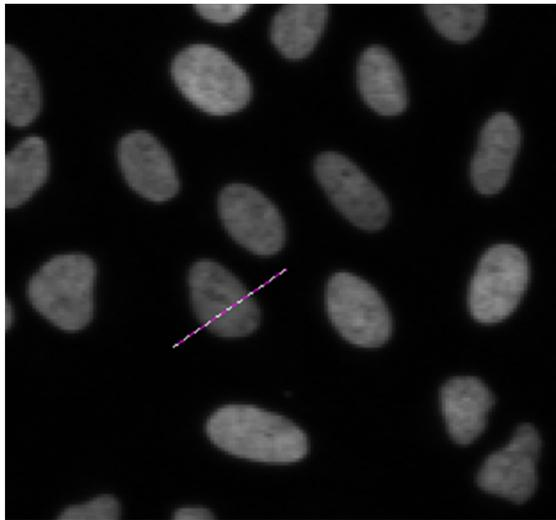
NOTE The above suggestions are only meant to guidelines for starting values. You will need to test and optimize these values.



Determining Intensity Above Local Background

The Intensity above local background is used for finding the cells:

- This value should be set to the intensity difference between a dim cell and its local background
- Using the **Region** Tools, draw a line region through a cell as shown below
- In the main menu, select **Measure > Intensities > Linescan**
- Measure the intensity change on the y-axis, then divide by 4
- Enter this result in the **Intensity above local background** box



Use the difference in intensity values to set **Intensity above local background**



Test Adaptive Acquisition Settings

Adaptive Acquisition

Minimum sites to visit: 2

Test Segmentation

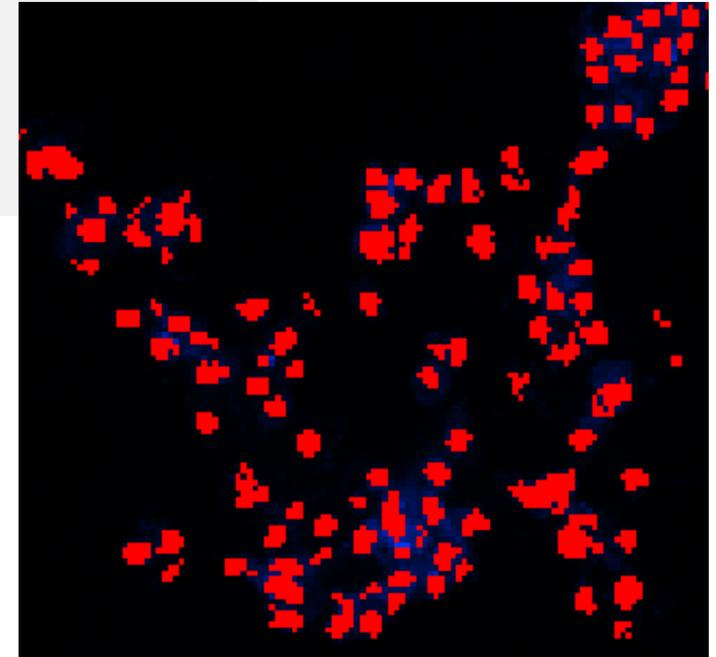
Wavelength: W1 - DAPI

Nuclei count: 43

Approximate width: 5 to 10 μm

Intensity above local background: 100 gray levels

Cell Count per well: 50

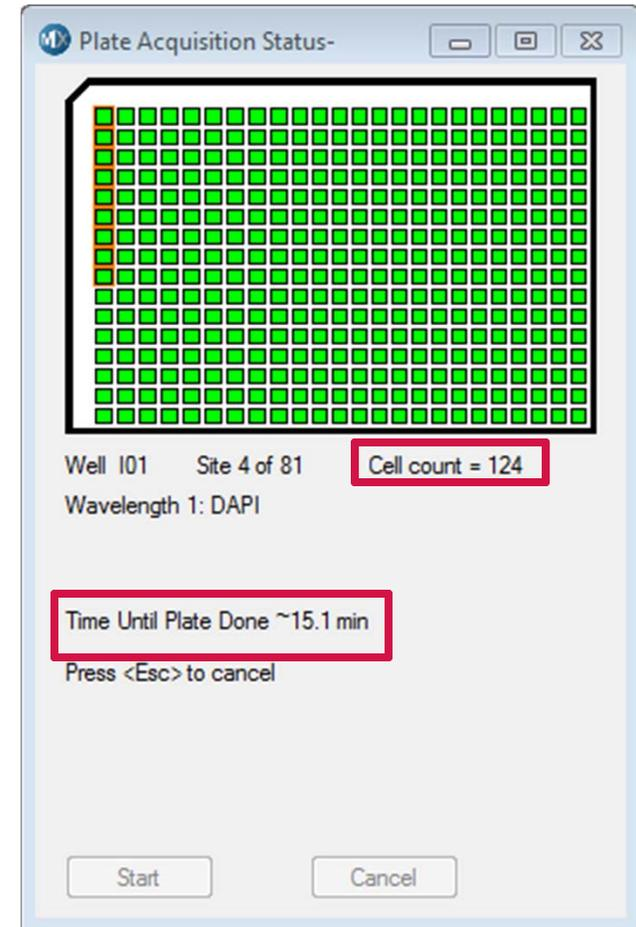


13. Click on the **Test Segmentation** button
 - An image of the selected wavelength will appear with a red segmentation overlay and the Nuclei (cell) count will be shown under the **Test Segmentation** button
14. If necessary, adjust the width and intensity settings to accurately find the cells
15. Continue to optimize settings for your plate as you would for normal acquisition and acquire your plate



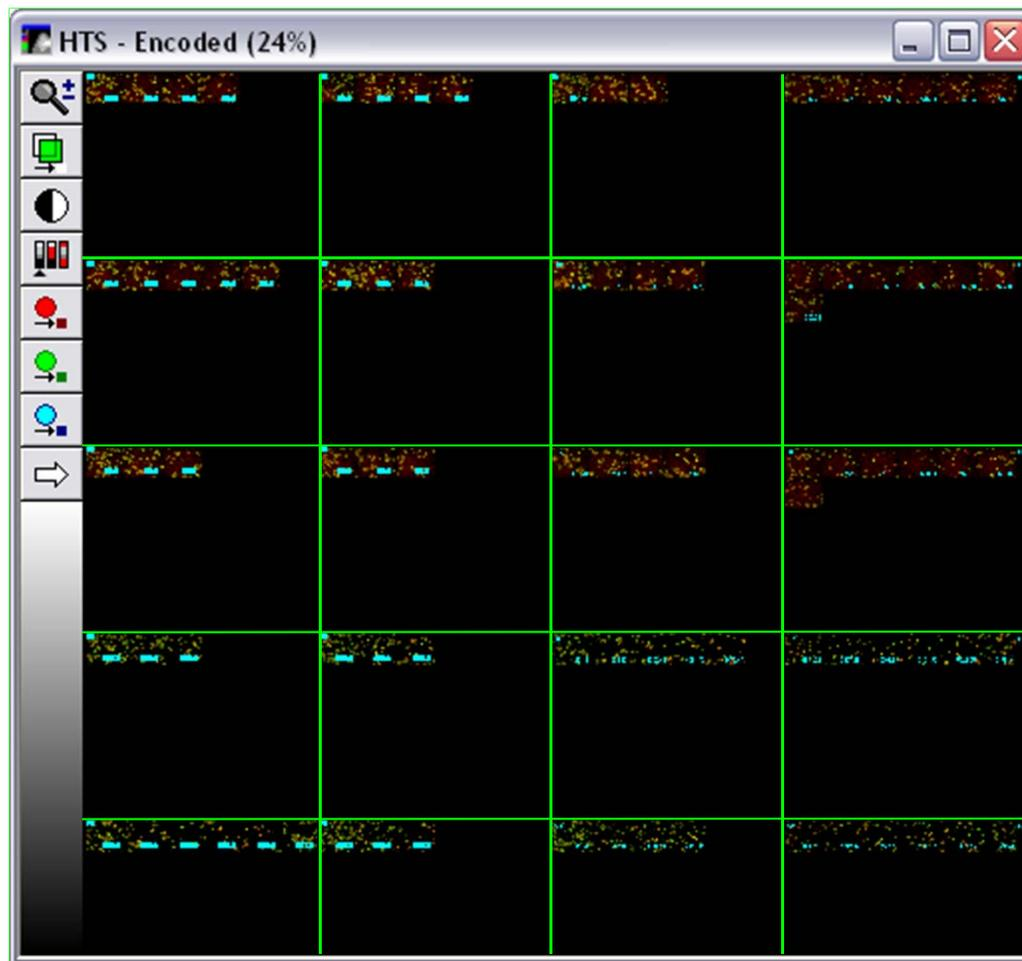
Acquiring with Adaptive Acquisition Enabled

- During plate acquisition, the Cell count will be shown on the **Plate Acquisition Status** dialog
- The **Time Until Plate Done** value should not be relied upon because it will vary considerably after each well is completed



Viewing Images in Review Plate Data

When you view thumbnail images through the **Review Plate Data** dialog, you will see that each well has a different number of sites acquired



Support Resources

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





ADVANCING PROTEIN AND CELL BIOLOGY