

MSD Assays

QUICKSTART guide for reading a plate
using MSD SECTOR[®] Imager

This package insert must be read in its entirety before using this product.

FOR RESEARCH USE ONLY.

NOT FOR USE IN DIAGNOSTIC OR THERAPEUTIC PROCEDURES.

Meso Scale Discovery

A division of Meso Scale Diagnostics, LLC.
9238 Gaither Road
Gaithersburg, MD 20877 USA
www.mesoscale.com

Meso Scale Discovery, Meso Scale Diagnostics, MSD, MSD (design), MULTI-ARRAY, MULTI-SPOT, SULFO-TAG and SECTOR are trademarks of Meso Scale Diagnostics, LLC.

© 2008 Meso Scale Discovery a division of Meso Scale Diagnostics, LLC. All rights reserved.



Reading a plate using MSD SECTOR Imager

Reading a plate using MSD SECTOR Imager



Check the Taskbar at the bottom of the screen for MSD Discovery Workbench®. If the program is not already opened, double-click the MSD Discovery Workbench icon on the desktop.



Click on the **SECTOR Imager icon** on the Toolbar.

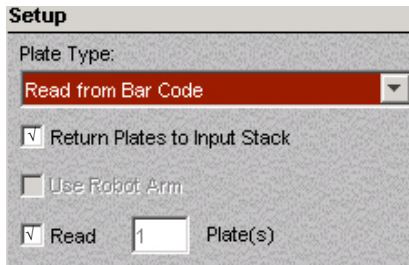
The SECTOR Imager splash screen will display while the instrument is initializing.

To read a single plate:

Ensure that the single plate adaptor is in the right-hand stack position
Load the MSD microplate into the single plate adaptor.

Note: Microplates can be loaded in either orientation. The software will automatically correct for differences in orientation provided that the barcode is intact.
Normal orientation is when well A1 is closest to the instrument (A1 at the top right hand corner).

To read a whole plate:



Select **“Read from Bar Code”** from the Plate Type menu.

Select the **“Return Plates to Input Stack”** checkbox.
Note: If left unchecked, the microplate will be transferred to the output stack on the left front side of the Sector Imager after it has been read.

Select the **Read Plate(s) checkbox** and enter **“1”** in the text field.

Click **Run** at the bottom of the SECTOR Imager window.

The Run Options window displays.

Verify the **“Export”** format choice and make changes if necessary (See User Manual for details).

Enter **RUN Name** – this is the file name under which the data will be saved.

Select **Output Path** – this is the location where the results file will automatically be saved.

Click **OK**. The plate read starts. It takes approximately 1minute (SI6000) or 3 minutes (SI2400) to read one plate.

You do not need to close this window to run another plate – simply make the appropriate selections as indicated above and proceed to read the next plate.

