

MSD Assays

Use and Storage of a Partial Plate

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.



Use and Storage of a Partial Plate

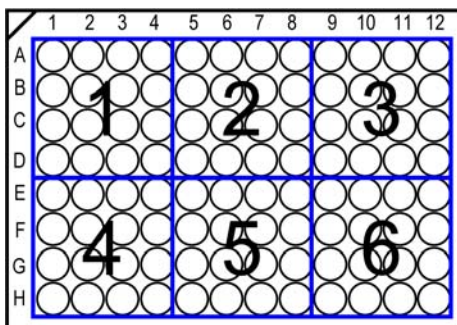
use and storage of a partial plate

Assays can be performed on parts of an MSD plate that can be then stored for later use. It is important to keep the unused wells dry during the assay. For proper use of partial plate the following recommendations should be followed:

1. The layout of the assay should be done within the boundary of each sector (see next page).
2. The amount of each component mixed should be adjusted for the fraction of the plate to be used.
3. The MSD plate packaging (bag and desiccant pouch) should be retained.
4. The unused sector(s) should be covered with an adhesive plate seal before starting the assay.
5. After addition of Read Buffer, the plate seal should be removed before the plate is placed into the SECTOR instrument.
6. To read a partial plate, the plate type should be chosen from the pull-down menu (the plate type is indicated on the packaging of the plate) in order for the partial plate layout to appear on the screen.
7. Select the Read Plate(s) checkbox and enter "1" in the text field.
8. Select the sector(s) in use and read the plate.
9. Following the partial plate read, cover the unused section(s) with an adhesive plate seal and remove the Read Buffer from the used wells. Tap out all residual fluid onto paper towels.
10. Return the sealed plate to the MSD plate bag (containing the desiccant pack) and place within a sealable plastic bag. Remove as much air from the bag as possible and seal.
11. Store plate for up to 2 weeks at 4°C.

Plate sector maps for SECTOR 6000 Imager and SECTOR 2400 Imager systems. Partial plate reads should use all wells with a specific sector and used sectors should be protected with an adhesive plate seal during the partial-plate assay.

SECTOR 6000 Imager Sector Map



SECTOR 2400 Imager Sector Map

