

How to Create a Kit Layout in DISCOVERY WORKBENCH® 4.0

Kit layouts are specific to a plate type (wells/plate and spots/well). You can create a new kit layout by modifying a pre-existing layout and saving it with a new name or by starting with a blank plate.

To Create a New Kit Layout:

- 1) Open the Kit Layout Editor (Tools → Kit Layout Editor).
- 2) Select the Create New Kit Layout icon (Figure 1). Give the layout a name, choose your plate type, and select a category for the assay/panel (Figure 2).
- 3) Right-click a spot to bring up the menu (Figure 3).
- 4) Select Assign Assay, then select the assay from the Select Assay window (Figure 4). The assay lists can be filtered or sorted. To enter an assay manually, select Create Assay (Figure 3).
- 5) Repeat for each spot.
- 6) Click the Save icon (Figure 1) to save the new kit layout on the Custom Layouts tab.

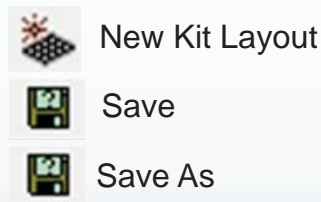


Figure 1: Kit Layout Editor icons

To Create a Kit Layout from an Existing Layout:

- 1) Open the Kit Layout Editor (Tools → Kit Layout Editor).
- 2) Select an existing catalog or custom kit layout that matches your plate configuration.
- 3) Right-click a spot to bring up the menu (Figure 3).
- 4) Select Assign Assay, then select the assay from the Select Assay window (Figure 4). The assay lists can be filtered or sorted. To enter an assay manually, select Create Assay (Figure 3).
- 5) Repeat for each spot.
- 6) Click the Save As icon (Figure 1) and type in a name to save the new kit layout. The new kit layout will be saved on the Custom Layouts tab.

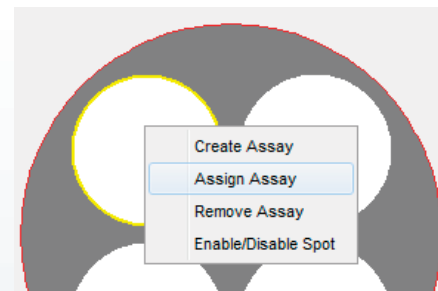


Figure 3: Assigning an assay to a spot

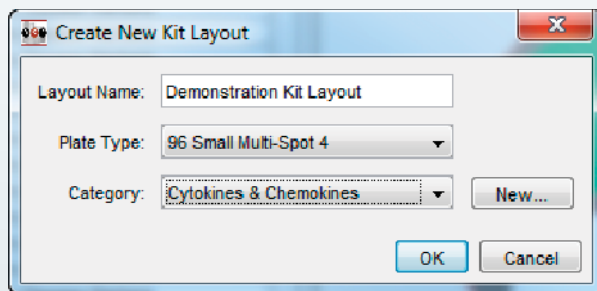


Figure 2: Create New Kit Layout window

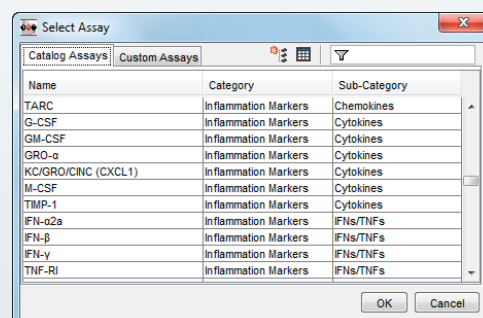


Figure 4: Select Assay window

©2013 Meso Scale Diagnostics, LLC. All Rights Reserved. MESO SCALE DISCOVERY, MESO SCALE DIAGNOSTICS, WWW.MESOSCALE.COM, MSD, MSD (DESIGN), DISCOVERY WORKBENCH, QUICKPLEX, MULTI-ARRAY, MULTI-SPOT, SULFO-TAG, SECTOR, SECTOR HTS, SECTOR PR, SMALL SPOT (DESIGN), 96 WELL 1, 4, 7, & 10-SPOT (DESIGNS), 384 WELL 1 & 4-SPOT (DESIGNS) and SPOT THE DIFFERENCE are trademarks and/or service marks of Meso Scale Diagnostics, LLC.