

Creating an Experiment in DISCOVERY WORKBENCH® 4.0

Date	Owner	Plate Type	Kit Layout	Long-side B... Code	Short-side ... Code	Manufacturer Bar Code	Quarantined	Imported	Creation Time	Hi
05/22/2012 1...	MSD Operator	96 Multi-Spot 7	Imported Lay...			29A3SAH038		true	05/22/2012 1...	
05/22/2012 1...	MSD Operator	96 Multi-Spot 7	Imported Lay...			29A3SAV082		true	05/22/2012 1...	
05/22/2012 1...	MSD Operator	96 Multi-Spot 7	Imported Lay...			29A3SAP113		true	05/22/2012 1...	
06/26/2012 0...	MSD Analyst	96 Multi-Spot	Human Bro I			2BCWZAK021			06/26/2012 0...	
06/26/2012 0...	MSD Analyst	96 Multi-S				A011024001			06/26/2012 0...	
06/26/2012 0...	MSD Analyst	96 Multi-S				B121019004			06/26/2012 0...	
06/26/2012 0...	MSD Analyst	96 Multi-S				25CWZAD021			06/26/2012 0...	
06/26/2012 0...	MSD Analyst	384				5101005005			06/26/2012 0...	

To create an experiment from a plate layout:

- 1) Launch DISCOVERY WORKBENCH software.
- 2) Go to the Plate Data History, and select the Plates (Table) or Plates (Tree) tab.
- 3) Right click on the plate you wish to analyze and select Analyze Plates.
- 4) You will be prompted to select a plate layout; only plate layouts of the correct well/spot configuration will be shown. Choose your plate layout and click OK.
- 5) This will create an experiment.

To create an experiment from a template:

- 1) Launch DISCOVERY WORKBENCH software.
- 2) Go to the Plate Data History, and select the Plates (Table) or Plates (Tree) tab.
- 3) Right click on the plate you wish to analyze and select Analyze Plates Using Template.
- 4) You will be prompted to select a Template; only templates of the correct well/spot configuration will be shown. Choose your template and click OK.
- 5) This will create an experiment using the information in the template, including the saved plate layout, table and plot formatting, and profiles.