

Releasing a Quarantined Plate in DISCOVERY WORKBENCH® 4.0

- 1) Launch DISCOVERY WORKBENCH software.
- 2) Open the Plate Data History and select the Plate (Tree) or Plate (Table) tab.
- 3) Locate a quarantined plate (Figures 1 and 2). Right click on the quarantined plate, and choose Release Plate.
- 4) The software will prompt you with what is required to release the plate.
 - a. Valid MSD data integrity key (if the bar code is expired or was unreadable/missing/already used).
 - b. Valid Kit Layout (if a kit layout for a different plate type was chosen).
- 5) Provide the requested information. The plate will open, and the software will prompt you to export the raw data, which is normally done automatically.
- 6) The run will be released and can be viewed, analyzed, and re-exported.

Note: A quarantined plate can also be released immediately after being read by providing a valid data integrity key or kit layout (Figure 3).

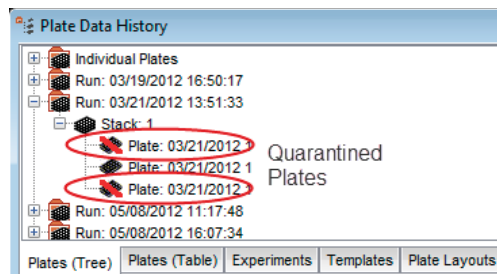


Figure 1. Plate Data History, Plate (Tree) view

Date	Owner	Plate Type	Kit Layout	Long-side B... Code	Short-side ... Code	Manufacturer Bar Code	Quarantined	Imported	Creation T
06/26/2012 0...	MSD Analyst	96 Multi-Spot...	Human Pro-L...			B121019004			06/26/2012
06/26/2012 0...	MSD Analyst	96 Multi-Spot 4	Human Pro-L...			25CWZAD021			06/26/2012
06/26/2012 0...	MSD Analyst	384	Standard			5101005005	Expired bar code		06/26/2012
06/26/2012 1...	MSD Analyst	96 Multi-Spot...	Standard			2BCWZAK021	Previously read plate		06/26/2012
06/26/2012 1...	MSD Analyst	96 Multi-Spot...	Standard			2BCWZAK021	Previously read plate		06/26/2012
06/26/2012 1...	MSD Analyst	96 Multi-Spot 7	Human Pro-L...			A01102400			06/26/2012
06/26/2012 1...	MSD Analyst	96 Multi-Spot 4	Standard			25CWZAD021	Invalid kit layout sele		06/26/2012
05/22/2012 1...	MSD Operator	96 Multi-Spot 7	Imported Lay...			29A3SAH038		true	05/22/2012
05/22/2012 1...	MSD Operator	96 Multi-Spot 7	Imported Lay...			29A3SAV082		true	05/22/2012
05/22/2012 1...	MSD Operator	96 Multi-Spot 7	Imported Lay...			29A3SAP113		true	05/22/2012

Figure 2. Plate Data History, Plate (Table) view

The plate has been read, but the data has been quarantined for certification. The data integrity records indicate that this plate has been previously used. Reusing a plate may generate compromised data.

If you would still like to release the data, you must enter a new data integrity key (barcode) in the box below. To obtain a key, please follow one of these options.

A. Go to <http://www.mesoscale.com/CatalogSystemWeb/WebRoot/products/ditkey.aspx?plate-type=96+Multi-Spot+10&serial=ST> and generate a new data integrity key. You will need to provide your email address.

B. Contact MSD Customer Service for assistance:
+1 301.947.2085 or CustomerService@mesoscale.com

Figure 3. Instructions for releasing a plate