



CERTIFICATE OF ANALYSIS

COA - 03676

Product Description: MULTI-ARRAY® 96-well High Bind Plate

Catalog #/Size: L15XB-1/ 1 plate; L15XB-3/ 10 plates; L15XB-6/ 100 plates; L15XB-7/ 500 plates

Lot Number: 22B12A

Expiration Date: 30 NOV 2013

Statement: This product was manufactured and tested according to approved MSD documents. This lot meets MSD's specifications.

BON (Electrical resistivity): < 200 microohms in all circuits

Leak test: No leaks

Free Tag 15,000 test

ECL (mean ± tolerance): 27,400 ± 5,480 cts

Max Average Intraplate CV for any spot: 3%

Maximum Intraplate CV for any plate: 7%

Max Interplate CV for any spot: 7%

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Preparer	Rolanda Spratley	Biologics/Chemistry		18 JUN 2013
Review/Approval	Michael Eager	Quality Control		05 JUL 2013