



## CERTIFICATE OF ANALYSIS

COA - 04014

**Product Description:** MULTI-ARRAY<sup>®</sup> 96-well High Bind Plate Pack

**Catalog #/Size:** L15XB-1 / 1 Plate; L15XB-3 / 10 Plates; L15XB-6 / 100 Plates; L15XB-7 / 500 Plates

**Lot Number:** 22D0WA

**Expiration Date:** 30 JUN 2015

**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  $\pm$  tolerance): 27 400  $\pm$  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
<b>Preparer</b>	Rolanda Spratley	Biologics/Chemistry		03 JAN 2014
<b>Review/Approval</b>	Michael Eager	Quality Control		06 JAN 2014