



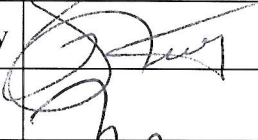

CERTIFICATE OF ANALYSIS

COA - 03463

**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:** 22C0PA  
**Expiration Date:** October 2014  
**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
<b>Preparer</b>	Rolanda Spratley	Biologics/Chemistry		15 Jan 13
<b>Review/Approval</b>	Michael Eager	Quality Control		15 JAN 2013