



CERTIFICATE OF ANALYSIS

COA – 06459

**Product Description:** MULTI-ARRAY® 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:** 22F0QA

**Expiration Date:** 30 NOV 2017

**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): 23 461 ± 4, 692 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.**

	<b>Name</b>	<b>Function</b>	<b>Signature</b>	<b>Date</b>
<b>Preparer</b>	Rolanda Spratley	Biologics/Chemistry		05 FEB 2016
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance		05 FEB 2016