



CERTIFICATE OF ANALYSIS

COA - 03489

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: 22C0NA
Expiration Date: August 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
Preparer	Rolanda Spratley	Biologics/Chemistry		07 FEB 13
Review/Approval	Michael Eager	Quality Control		12 FEB 2013