



CERTIFICATE OF ANALYSIS

COA - 03828

Product Description: MULTI-ARRAY[®] 96-well Plate Pack

Catalog #/Size: L15XA-1 / 1 Plate; L15XA-3 / 10 Plates; L15XA-6 / 100 Plates; L15XA-7 / 500 Plates

Lot Number: 21D1DA

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

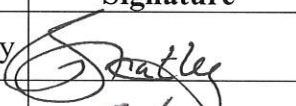

ECL (mean \pm tolerance): 15 000 \pm 3,000 cts

Max Average Intraplate CV for any spot: 4%

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		08 AUG 2013
Review/Approval	Michael Eager	Quality Control		26 AUG 2013