



CERTIFICATE OF ANALYSIS

COA - 06149

Product Description: MULTI-ARRAY[®] 96, PR,1 Spot Std,Plate Assy, Lipless Top

Catalog #/Size: 3-2010-120002-F

Lot Number: 01E01A

Expiration Date: 30 APR 2016

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: 3%

Maximum Intraplate CV for any plate: 7%

Max Interplate CV for any spot: 7%

96-1 LS Spot STD

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Preparer	Rolanda Spratley	Biologics/Chemistry		28 OCT 2015
Review/Approval	Matthew Chorley	Quality Assurance		29 OCT 2015