



CERTIFICATE OF ANALYSIS

COA - 05435

Product Description: MULTI-ARRAY[®] 96-well 5m Spot High Bind Plate Pack

Catalog #/Size: L45XB-1 / 1 Plate; L45XB-3 / 10 Plates; L45XB-6 / 100 Plates; L45XB-7 / 500 Plates

Lot Number: 24E07A

Expiration Date: 30 JUN 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): 27040 \pm 5408 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		19 MAY 2015
Review/Approval	Anastasia Belilovets	Quality Assurance		20 MAY 2015