



CERTIFICATE OF ANALYSIS

COA - 05411

Product Description: MULTI-ARRAY® 96-well, 5mm Spot Plate Pack

Catalog #/Size: L45XA-1 / 1 Plate; L45XA-3 / 10 Plates; L45XA-6 / 100 Plates; L45XA-7 / 500 Plates

Lot Number: 23E09A

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 ± tolerance): 15 000 ± 3,000 cts

Max Average Intraplate CV for any spot: 3%

Maximum Intraplate CV for any plate: 6%

Max Interplate CV for any spot: 6%

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

	Name	Function	Signature	Date
Preparer	Rolanda Spratley	Biologics/Chemistry		24 APR 2015
Review/Approval	Anastasia Belilovets	Quality Assurance		24 APR 2015