



CERTIFICATE OF ANALYSIS

COA - 05048

Product Description: MULTI-ARRAY[®] MA6000 384 HB Plate Pack

Catalog #/Size: L21XB-1 / 1 Plate; L21XB-4 / 15 Plates; L21XB-6 / 120 Plates; L21XB-7 / 510 Plates

Lot Number: 06D0FA

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): N/A

Max Average Intraplate CV for any spot: 5%

Maximum Intraplate CV for any plate: 7%

Max Interplate CV for any spot: 8%

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

	Name	Function	Signature	Date
Preparer	Rolanda Spratley	Biologics/Chemistry		13 JAN 2015
Review/Approval	Anastasia Belilovets	Quality Control		19 JAN 2015