



CERTIFICATE OF ANALYSIS

COA - 04098

Product Description: MA6000 384 Plate Pack

Catalog #/Size: L21XA-1 / 1 Plate; L21XA-4 / 15 Plates; L21XA-6 / 120 Plates; L21XA-7 / 510 Plates

Lot Number: 05D04A

Expiration Date: 31 MAY 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): 18 000 \pm 3,600 cts

Max Average Intraplate CV for any spot: 5%

Maximum Intraplate CV for any plate: 9%

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		24 FEB 2014
Review/Approval	Michael Eager	Quality Control		24FEB2014