



## CERTIFICATE OF ANALYSIS

COA - 04102

**Product Description:** MA2400 384 Plate Pack

**Catalog #/Size:** L25XA-1 / 1 Plate; L25XA-4 / 15 Plates; L25XA-6 / 120 Plates; L25XA-7 / 510 Plates

**Lot Number:** 11D01A

**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): 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
<b>Preparer</b>	Rolanda Spratley	Biologics/Chemistry		28 JAN 2014
<b>Review/Approval</b>	Michael Eager	Quality Control		10 FEB 2014