



CERTIFICATE OF ANALYSIS

COA - 03861

Product Description: MULTI-ARRAY® PR HB Plate Pack
Catalog #/Size: L13XB-1 / 1 Plate; L13XB-3 / 10 Plates; L13XB-6 / 100 Plates; L13XB-7 / 500 Plates
Lot Number: 02D01A
Expiration Date: 30 APR 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 ± tolerance): 30 000 ± 6,000 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		13SEP2013
Review/Approval	Michael Eager	Quality Control		17SEP2013