



CERTIFICATE OF ANALYSIS

COA - 03488

Product Description: MULTI-ARRAY® 96-well, 5mm Spot Std Plate Pack
Catalog #/Size: L45XA-1 / 1 Plate; L45XA-3 / 10 Plates; L45XA-6 / 100 Plates; L45XA-7 / 500 Plates
Lot Number: 23C0XA
Expiration Date: October 2014
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.

Table with 5 columns: Name, Function, Signature, Date. Rows include Preparer (Rolanda Spratley) and Review/Approval (Michael Eager).