



CERTIFICATE OF ANALYSIS

COA - 03870

Product Description: MULTI-ARRAY® 96-well High Bind Plate Pack
Catalog #/Size: L15XB-1 / 1 Plate; L15XB-3 / 10 Plates; L15XB-6 / 100 Plates; L15XB-7 / 500 Plates
Lot Number: 22D0GA
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 ± tolerance): 27 400 ± 5,480 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.

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