



CERTIFICATE OF ANALYSIS

COA – 07088

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: 22G0HA

Expiration Date: 31 MAY 2018

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): 23 461 \pm 4, 692 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 | | 07 JUL 2016 |
| Review/Approval | Dawn Stollar | Quality Assurance | | 07 JUL 2016 |