



CERTIFICATE OF ANALYSIS

COA - 06641

Product Description: MULTI-SPOT® 96, SI, 4 Spot Std Lipless Top Plate

Catalog #/Size: 3-2024-120002 / 1 Plate

Lot Number: 25F3YA

Expiration Date: 30 JUN 2017

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: 5%

Maximum Intraplate CV for any plate: 7%

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 |
|------------------------|------------------|---------------------|---|-------------|
| Preparer | Rolanda Spratley | Biologics/Chemistry |  | 15 MAR 2016 |
| Review/Approval | Dawn Stollar | Quality Assurance |  | 15 MAR 2016 |