



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: U-PLEX[®] 10-Assay, 96-Well SECTOR[®] Plate

Catalog Number: N05235A-1

Lot Number: Z00U0040

Expiration Date: 31 DEC 2019

Storage Condition: 2-8°C

Acceptance Data:

| Metric | Specifications | All Spots |
|----------------------------------|----------------|-----------|
| Average Intraplate CV | ≤ 10.0% | Pass |
| Average Interplate CV | ≤ 13.0% | Pass |
| Signal Range at Binding Capacity | ≤ 30.0% | Pass |

Statement: The above plate lot is intended for Research Use Only. Not for use in Diagnostic Procedures.

| | Name | Function | Signature | Date |
|------------------------|--------------|-------------------|---------------------|-------------|
| Review/Approval | Dawn Stollar | Quality Assurance | <i>Dawn Stollar</i> | 13 SEP 2017 |