



CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-07229

Product Description: U-PLEX[®] 2-Assay, 96-Well SECTOR[®] Plate
Catalog Number: N05227A-1
Lot Number: Z00U0023A
Expiration Date: 31 JAN 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 was tested using MSD procedures and conforms to MSD's specifications. It is intended for Research Use Only. Not for use in Diagnostic Procedures.

| | Name | Function | Signature | Date |
|-----------------|------------------------|----------------------|----------------------------|-----------|
| Review/Approval | Vanessa Castleberry | Quality Assurance | <i>Vanessa Castleberry</i> | 19AUG2016 |