



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** U-PLEX® 6-Assay, 96-Well SECTOR® Plate

**Catalog Number:** N05231A-1

**Lot Number:** Z00U0047E

**Expiration Date:** 31 MAR 2020

**Storage Condition:** 2-8°C

### Acceptance Data:

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

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

|                        | Name         | Function          | Signature           | Date        |
|------------------------|--------------|-------------------|---------------------|-------------|
| <b>Review/Approval</b> | Dawn Stollar | Quality Assurance | <i>Dawn Stollar</i> | 01 DEC 2017 |