



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** U-PLEX<sup>®</sup> 7-Assay, 96-Well SECTOR<sup>®</sup> Plate

**Catalog Number:** N05232A-1

**Lot Number:** Z00U0057F

**Expiration Date:** 31 AUG 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        |
|-----------------|--------------|-------------------|---------------------|-------------|
| Review/Approval | Dawn Stollar | Quality Assurance | <i>Dawn Stollar</i> | 20 MAR 2018 |