



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: U-PLEX[®] 3-Assay, 96-well SECTOR[®] Plate

Catalog Number: N05228A-1

Lot Number: Z00U0141B

Expiration Date: 31 AUG 2023

Storage Temperature: 2-8°C

Acceptance Data:

| Metric | Specifications | All Spots |
|---------------------------|----------------|-----------|
| Average Intraplate CV | $\leq 10.0\%$ | Pass |
| CV of Intraplate Averages | $\leq 13.0\%$ | Pass |

This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

| | Name | Function | Signature | Date |
|-----------------|------------------|----------|-----------|-------------|
| Review/Approval | Xiomara Talavera | Quality | | 25 JUN 2021 |