



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** U-PLEX<sup>®</sup>10-Assay, 96-well SECTOR<sup>®</sup> Plate

**Catalog Number:** N05235A-1

**Lot Number:** Z00U0139


**Expiration Date:** 31 AUG 2023

**Storage Temperature:** 2-8°C

### Acceptance Data:

Metric	Specifications	All Spots
Average Intraplate CV	≤ 10.0%	Pass
CV of Intraplate Averages	≤ 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		08 OCT 2021