



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** U-PLEX<sup>®</sup> 3-Assay, 96-well SECTOR<sup>®</sup> Plate

**Catalog Number:** N05228A-1

**Lot Number:** Z00U0128B


**Expiration Date:** 31 MAR 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	Karen Hamilla	Quality		23 FEB 2021