



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** U-PLEX® 2-Assay, 96-Well SECTOR® Plate

**Catalog Number:** N05227A-1

**Lot Number:** Z00U0053A

**Expiration Date:** 30 JUN 2020

**Storage Condition:** 2-8°C

### Acceptance Data:

Metric	Specifications	All Spots
Average Intraplate CV	≤ 10.0%	Pass
Average Interplate CV	≤ 13.0%	Pass
Signal Range at Binding Capacity	≤ 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>	30 JAN 2018