



### CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** U-PLEX® 10-Assay, 96-Well SECTOR® Plate

**Catalog Number:** N05235A-1

**Lot Number:** Z00U0033

**Expiration Date:** 31 OCT 2019

**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
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	05 JUN 2017