



## 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:** Z00U0076

**Expiration Date:** 30 APR 2021

**Storage Condition:** 2-8°C

### Acceptance Data:

Metric	Specifications	All Spots
Average Intraplate CV	$\leq 10.0\%$	Pass
Average Interplate CV	$\leq 13.0\%$	Pass
Signal Range at Binding Capacity	$\leq 30.0\%$	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>	04 JAN 2019