



## CERTIFICATE OF ANALYSIS

Page 1 of 1

COA-07423

**Product Description:** U-PLEX<sup>®</sup> 2-Assay, 96-Well SECTOR<sup>®</sup> Plate

**Catalog Number:** N05227A-1

**Lot Number:** Z00U0025A

**Expiration Date:** 31 MAR 2019

**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 was tested using MSD procedures and conforms to MSD's specifications. It 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>	19 OCT 2016