



## RFE MEMO


Page 1 of 1

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

**Catalog Number(s) / Size(s):** N05227A-1

**Lot Number:** Z00U0072A

Based on additional testing, the original date of expiration for this plate has been extended by 12 months. The expiration date of Z00U0072A has been extended to 28 FEB 2022.

Function	Name	Signature	Date
Quality	Karen Hamilla		14 MAY 2021

This memo supersedes all other statements regarding date of expiration for this product.



## CERTIFICATE OF ANALYSIS

Page 1 of 1

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

**Catalog Number:** N05227A-1

**Lot Number:** Z00U0072A

**Expiration Date:** 28 FEB 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
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	05 NOV 2018