



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: U-PLEX[®] 3-Assay, 96-Well SECTOR[®] Plate

Catalog Number: N05228A-1

Lot Number: Z00U0068B

Expiration Date: 31 JAN 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>	11 SEP 2018