



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: U-PLEX[®] 5-Assay, 96-Well SECTOR[®] Plate

Catalog Number: N05230A-1

Lot Number: Z00U0065D

Expiration Date: 31 DEC 2020

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
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	03 AUG 2018