



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: U-PLEX[®] 6-Assay, 96-well SECTOR[®] Plate

Catalog Number: N05231A-1

Lot Number: Z00U0146E


Expiration Date: 30 NOV 2023

Storage Temperature: 2-8°C

Acceptance Data:

Metric	Specifications	All Spots
Average Intraplate CV	≤ 10.0%	Pass
CV of Intraplate Averages	≤ 13.0%	Pass

This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Xiomara Talavera	Quality		08 OCT 2021