



CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog Number: N05228A-1

Lot Number: Z00U0175B

Expiration Date: 31 JUL 2024

Storage Temperature: 2-8°C

Acceptance Data:

Metric	Specifications	All Spots
Average Intraplate CV	$\leq 10.0\%$	Pass
CV of Intraplate Averages	$\leq 13.0\%$	Pass

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

	Name	Function	Signature	Date
Review/Approval	Xiomara Talavera	Quality		28 JUN 2022