



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: U-PLEX® 4-Assay, 96-Well SECTOR® Plate

Catalog Number: N05229A-1

Lot Number: Z00U0083C

Expiration Date: 31 AUG 2021

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	Dawn Stollar	Quality		24 APR 2019