



CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog Number: N05229A-1

Lot Number: Z00U0130C


Expiration Date: 30 APR 2023

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	Nithya Santhanam	Quality		24 Feb 2023